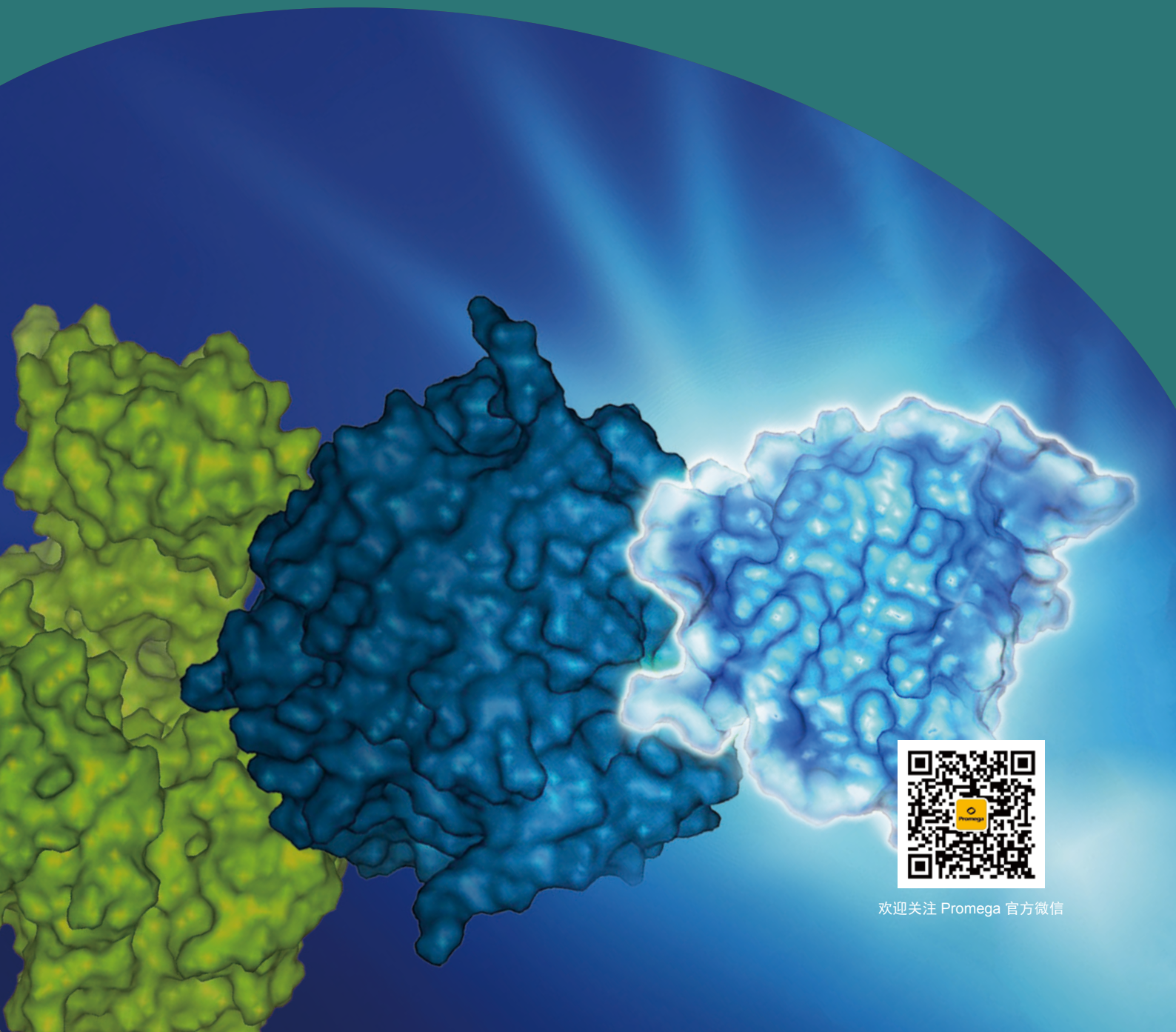


Dual Luciferase Reporter Technology

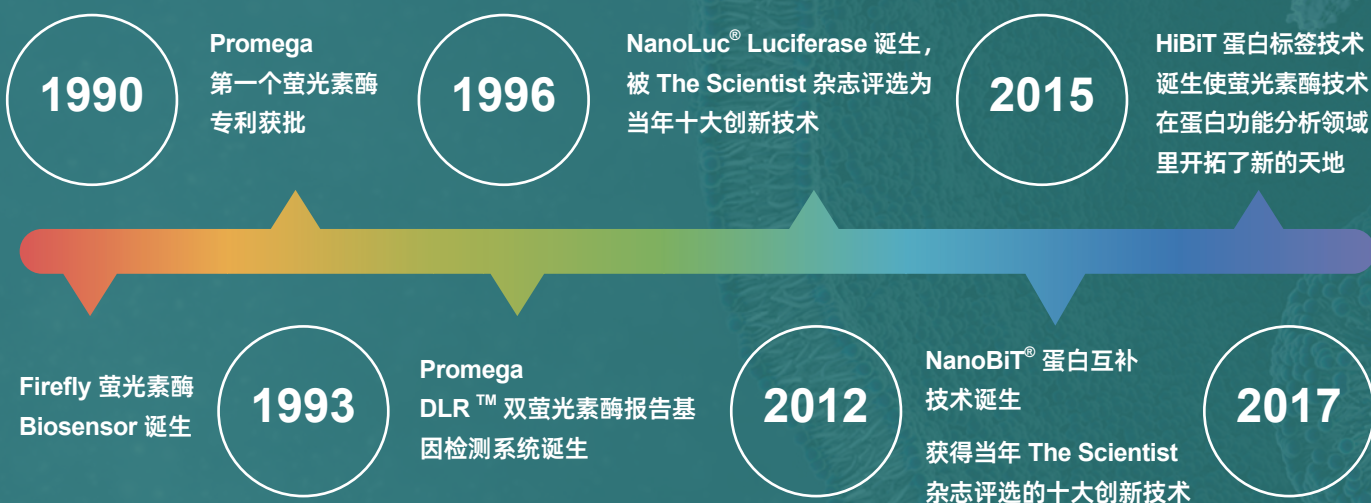
双萤光素酶报告基因解决方案



欢迎关注 Promega 官方微信

Promega 作为报告基因检测方案的最资深供应商，一直专注于基于生物发光的萤光素酶报告基因技术的开发。Promega 将一直致力于技术创新，为客户提供更为精良的研发工具和更完善的服务。

Promega 萤光素酶报告基因技术发展的时间轴



目 录

- 萤光素酶报告基因技术简介 3
- 双萤光素酶报告基因技术简介 4
- 双萤光素酶报告基因技术流程 5
- 双萤光素酶报告基因载体 6
- 双萤光素酶报告基因载体列表 8
- 双萤光素酶报告基因检测试剂 11
- 双萤光素酶报告基因检测相关产品..... 13
- 萤光素酶报告基因技术应用文献 15
- 产品列表 17

Promega 萤光素酶报告基因技术简介

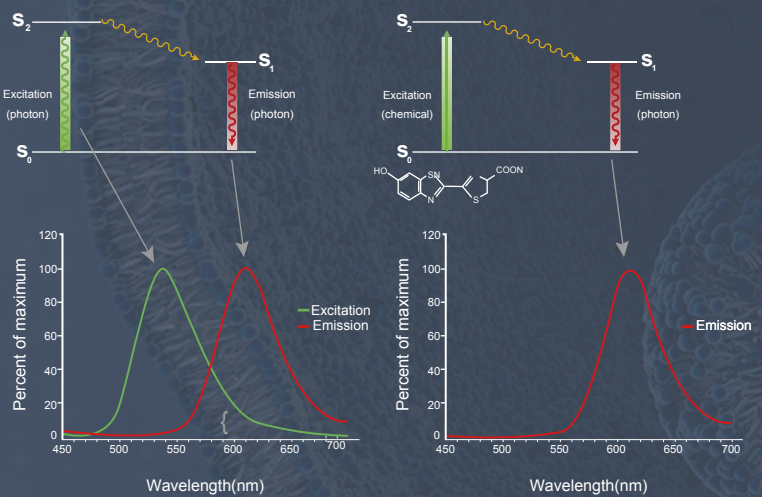
不同于荧光，生物发光依靠酶促反应产生自发光信号，无需激发，背景更低，灵敏度更高，不受化合物自发荧光干扰，更加适合作为细胞学分析手段。

目前除了广为人知的萤火虫萤光素酶报告基因，海肾萤光素酶报告基因以外，Promega 还最新推出了革新性的 NanoLuc[®] 报告基因，具有蛋白更小，表达更高，光信号更强等特点。

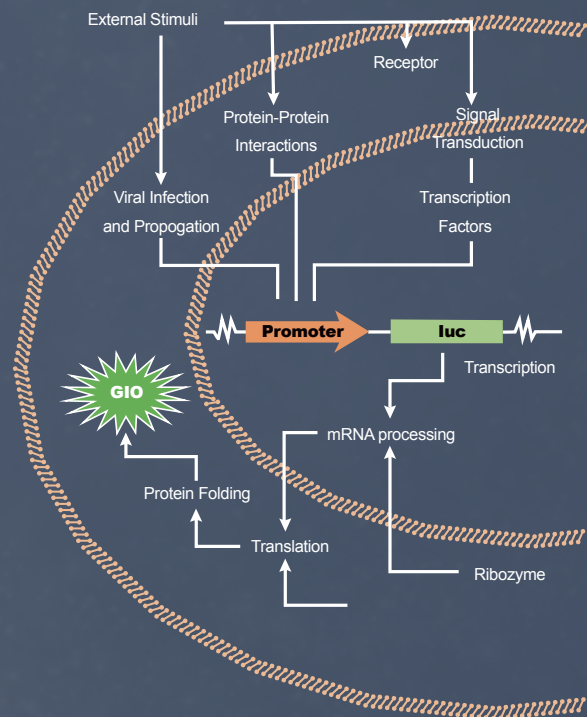
Promega 将一直致力于技术创新，为客户提供更为精良的研发工具和更完善的服务。

• 萤光素酶生物发光的特点：

- ✓ 无需激发：基于萤光素酶与底物的酶学反应产生自发光，无需激发光
- ✓ 背景低，灵敏度高：无激发光造成的背景，灵敏度更高，比荧光灵敏度高 100 倍
- ✓ 不受化合物自发荧光干扰：待测化合物经常会产生自发荧光，从而对基于荧光法的检测造成干扰，生物发光法则避免了这个问题



• 报告基因检测方案的应用



- ✓ 调控元件功能研究 (启动子, 增强子等)
- ✓ 细胞信号通路分析
- ✓ 膜受体和细胞内受体激活 / 结合研究
- ✓ 核受体研究
- ✓ siRNA/miRNA 研究
- ✓ 蛋白 - 蛋白相互作用
- ✓ 转录因子结果分析
- ✓ 病毒感染
- ✓ 干细胞分化
- ✓ 小动物活体成像
- ✓ 植物研究

双萤光素酶报告基因技术简介

• 什么是双萤光素酶报告基因检测？

是基于双萤光素酶报告基因载体构建、转染、表达，最终以检测系统来检测数据的一套研究方案，适用于多种应用方向。

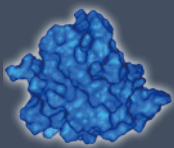
通常萤火虫 (Firefly) 萤光素酶报告基因载体作为目的片段克隆的载体，即主报告基因，海肾 (*Renilla*) 萤光素酶报告基因载体 (带组成型启动子) 作为内参载体。两个报告基因载体共转染，“内参对照”报告基因，用于归一化实验结果。对照报告基因可以控制细胞数量，细胞状态及转染效率造成的误差。

• 什么时候使用双报告基因检测？

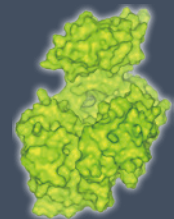
满足以下任意条件将需要进行双报告基因转染及检测，若均为否定答案则只需单报告基因转染及检测

- ✓ 需进行瞬时转染
- ✓ 需使报告基因信号减弱，比如研究负调控对刺激产生的细胞反应
- ✓ 需要共报告基因 (co-reporter) 以建立内对照或实现多重信号
- ✓ 实现最高的检测精度 (板内或板间误差更小)

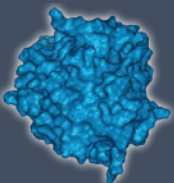
萤光素酶报告基因种类



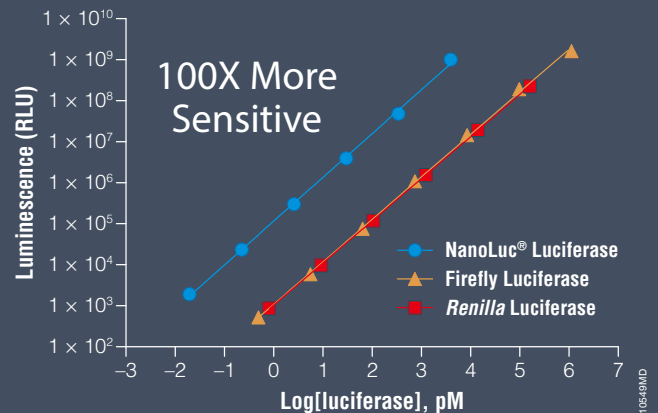
- NanoLuc® 萤光素酶 – NanoLuc® Luciferase
Promega 最新推出的专利报告基因，是一个经过基因工程改造的小分子酶 (19.1kDa)，配合革新型底物——Furimazine，可产生高强度、辉光型发光信号。生物发光反应不依赖 ATP，自发光背景低，光信号更强，同时抑制背景发光以获得最高检测灵敏度。既可用于主报告基因，也可用作内参报告基因。



- 萤火虫萤光素酶 – Firefly Luciferase
最经典常用的 Promega 生物发光报告基因，其产生的单体萤光素酶蛋白无需翻译后修饰，稳定性和表达效率更高。结合特别的快速应答技术，使其更适用于报告基因的应用，主要作为主报告基因。



- 海肾萤光素酶 – *Renilla* Luciferase
经典报告基因，来源于海洋中腔肠动物海肾，主要用于双报告基因检测中的内参对照报告基因。

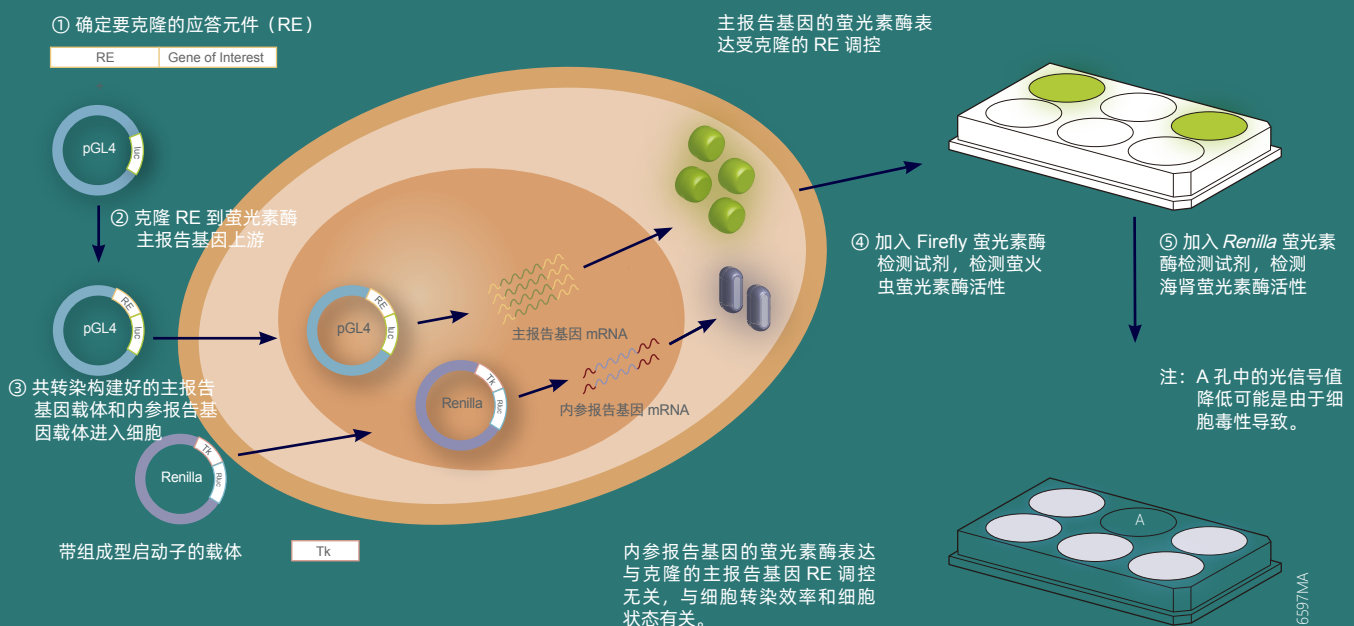


报告基因	NanoLuc®	Firefly	<i>Renilla</i>
来源	细角刺虾萤光素酶突变体	鞘翅目萤火虫	腔肠动物海肾
分子量	19kD	61kD	36kD
蛋白表达	细胞内或分泌型	细胞内	细胞内
颜色	蓝色	黄绿色	蓝色
底物	Furimazine	Luciferin	coelenterazine
灵敏度	+++++	++++	++++
线性范围	10 ¹⁰	10 ⁸	10 ⁸
主要应用	萤光素酶应用，融合蛋白表达，BRET，细胞信号通路分析	经典萤光素酶应用，细胞信号通路分析，调控元件功能型研究等	与 Firefly 共同使用做为内参报告基因

双萤光素酶报告基因技术流程

1 报告基因载体克隆构建	2 质粒提取, 细胞转染, 孵育	3 检测信号
<p>几种报告基因的的搭配选择: 主报告基因 / 内参报告基因</p> <ul style="list-style-type: none"> ④ NanoLuc®/Firefly ④ Firefly/NanoLuc® ④ Firefly/<i>Renilla</i> <p>Promega 报告基因载体工具:</p> <ul style="list-style-type: none"> ④ 骨架载体: 不带有任何应答元件, 有细胞内表达和分泌型表达可以选。 ④ 内参载体: 带有组成型启动子 ④ 预构建载体: 已预构建所需调控元件的载体 	<p>选择去除内毒素的质粒提取试剂盒</p> <ul style="list-style-type: none"> ④ 内毒素会影响细胞转染效果 <p>根据实验目的选择转染方法和转染试剂</p> <ul style="list-style-type: none"> ④ 转染方法: 瞬时转染或稳定转染 ④ 转染试剂: 转染效率高, 毒性低 ④ 稳转细胞: 已稳定转染所需目的基因或报告基因的细胞系 	<p>根据选择的报告基因选择检测试剂</p> <ul style="list-style-type: none"> ④ Firefly 与 NanoLuc® 组合: <ul style="list-style-type: none"> • NanoDLR™ ④ Firefly 与 <i>Renilla</i> 组合: <ul style="list-style-type: none"> • 非均质法 (DLR) • 均质法 (Dual-Glo™) <p>选择发光检测仪器</p> <ul style="list-style-type: none"> • 如选择非均质法检测试剂 (DLR) 通量高时就必须选择带有自动进样器的仪器。 • 选择高灵敏度检测仪。
<ul style="list-style-type: none"> • pGL4 系列萤光素酶载体 • pNL 系列萤光素酶载体 	<ul style="list-style-type: none"> • PureYield™ 质粒小提试剂盒 • ViaFect™ 转染试剂 • 商品化稳转细胞系 	<ul style="list-style-type: none"> • 检测试剂: 双报告基因检测试剂 (DLR, Dual-Glo™, NanoDLR™) • 检测仪器: GloMax® 系列检测仪

下图举例说明技术流程: 以 Firefly 萤光素酶报告基因 (Luc 或 Luc 2) 为主报告基因, 以 *Renilla* 萤光素酶报告基因 (RLuc) 为内参报告基因



双萤光素酶报告基因载体

Promega 拥有萤光素酶领域的专利技术，对萤光素酶工具的研究和开发处于全球领先的地位。pGL4 载体是继 pGL3 载体之后最新一代的载体。提供 NanoLuc[®]、Firefly 和 *Renilla* 三种萤光素酶报告基因载体。最新的 NanoLuc[®] 萤光素酶 pNL 系列载体，也是以 pGL4 载体为骨架设计，pNL 载体系列提供非融合和分泌型 Nluc 对照质粒。除骨架载体外，Promega 还提供预构建好各种信号通路启动元件的即用型载体。

• Promega 提供的报告基因载体种类：

<p>无启动子 / 含 minP 的报告基因载体 1</p> <p>用于转录水平的分析 (例如：启动子，启动子片段，信号转导研究)，NanoLuc[®] 系列有细胞内表达和分泌型两种可选。</p>	<p>应答元件 预构建载体 2</p> <p>近百种预构建好各种信号通路启动元件的即用型载体。可直接选购，无需自行构建。</p>	<p>报告基因 融合载体 3</p> <p>用于产生转录融合以研究转录 / 翻译后调控。用于感兴趣的蛋白和报告基因之间进行融合表达。</p>	<p>对照 内参载体 4</p> <p>选择带有组成型表达启动子的萤光素酶载体来作为内参。</p>
--	---	---	--

• 基因表达研究的报告基因载体设计：

转录调控 - 如：检测启动子活性



转录后调控 - 如：剪接 (Splicing)



转录后 / 翻译调控 - 如：检测 RNA 干扰 (RNAi)



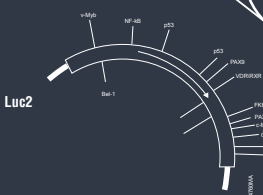
翻译后调控 - 如：检测蛋白稳定性



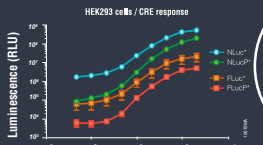
pGL4/pNL 载体主要特点

- 去除过氧化物酶靶标位点和隐秘的调控序列 - 减少非特异性表达。

载体骨架 更加优化



精确快速的 报告基因 应答反应

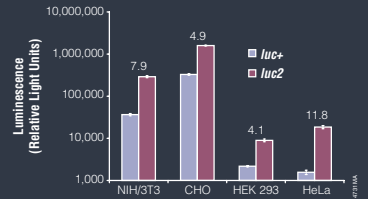


- 更加精确快速的报告基因应答反应 - pGL4/pNL 系列提供快速应答功能的载体，报告基因应答更快，生物相关性更好；带有快速应答功能的载体在报告基因缩写后会标有 P 或 CP。

Vector	Gene	Design
pGL4.10[luc2]	luc2	
pGL4.11[luc2P]	luc2P	
pGL4.12[luc2CP]	luc2CP	
pGL4.70[hRluc]	hRluc	

表达水平 更高 背景更低

- 系统性优化编码密码子 - 报告基因表达水平更高
- 背景更低，更灵敏 - 数据更出色



工具选择 更多

- 载体的选择 - 见报告基因载体列表。pGL4/pNL 载体的名称中均明确标示了其带有的主要元件。
- 商品化预构建信号通路调控元件载体 - 提供近百种萤光素酶报告基因和预构建元件骨架载体，减少自行构建的时间。

选择一个合适的载体组合

在 NanoLuc® 萤光素酶诞生以前，经典的萤光素酶双报告基因载体组合就是 Firefly 报告基因作为主报告基因，带组成型启动子的 *Renilla* 报告基因作为内参报告基因，NanoLuc® 萤光素酶报告基因诞生后，增加了这种组合的可选择性：

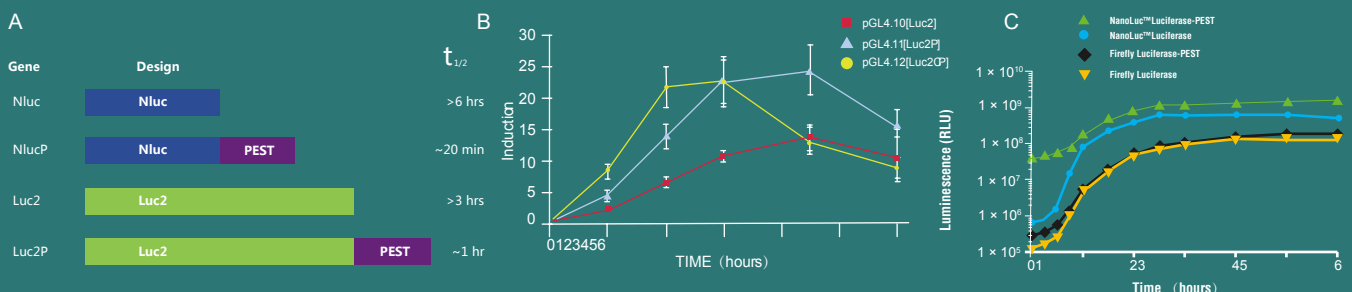
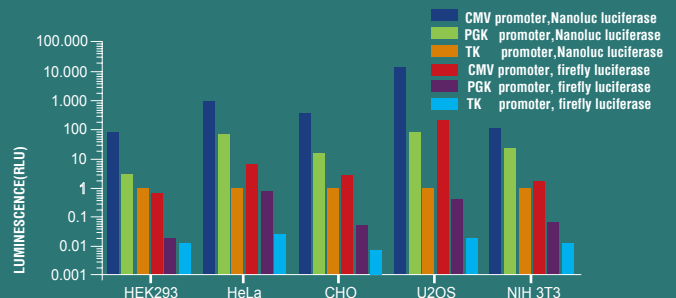
注意：Renilla 和 NanoLuc® 不能作为双报告基因组合使用。

- 主报告基因载体的选择：如果是基因调控研究，需要选择不带启动子或其他调控元件的骨架载体，根据组合不同，可选择 Firefly, NanoLuc® 中的任意一种。
- 内参报告基因载体的选择：带有表达组成型启动子的萤光素酶表达载体来作为内参，根据组合不同，内参报告基因可以选择 Firefly, NanoLuc® 或 *Renilla* 任意一种。推荐 pGL4 和 pNL 中带有不同强度的组成型启动子载体。

组合	主报告基因(含推荐载体举例)	内参报告基因(含推荐载体举例)	检测试剂
Firefly/ <i>Renilla</i>	Firefly pGL4.10[luc2] Vector (Cat.# E6651)	<i>Renilla</i> pGL4.74[hRluc/TK] Vector (Cat.# E6921) pGL4.73[hRluc/SV40] Vector (Cat.# E6911) pGL4.75[hRluc/CMV] Vector (Cat.# E6931)	DLR™ (Cat.# E1910系列) Dual-Glo® (Cat.# E2920系列)
NanoLuc®/Firefly	NanoLuc® pNL1.1[Nluc] Vector (Cat.# N1001)	Firefly pGL4.53[luc2/PGK] Vector (Cat.# E5011) pGL4.54[luc2/TK] Vector (Cat.# E5061) pGL4.13[luc2/SV40] Vector (Cat.# E6681)	NanoDLR™ (Cat.# N1610系列)
Firefly/NanoLuc®	Firefly pGL4.10[luc2] Vector (Cat.# E6651)	NanoLuc® pNL1.1.PGK[Nluc/PGK] Vector (Cat.# N1441) pNL1.1.TK[Nluc/TK] Vector (Cat.# N1501) pNL1.1.CMV[Nluc/CMV] Vector (Cat.# N1091)	NanoDLR™ (Cat.# N1610系列)

实验 Tips:

- 如果组合中使用 Nluc，则可以使用更高的报告基因载体用量比例，可能最高可以达到 1000:1 (Fluc:Nluc)。
- 同一代的内参载体之间，最主要的区别是携带的启动子不同，包括 CMV、SV40、TK 等。通常，内参基因表达的活性应当显著的高于背景组，同时，尽量小，确保不干扰主报告基因的表达。所以，我们建议选择内参载体时，首先考虑活性较弱的 TK 启动子驱动的载体，启动子强弱见图。如果通过文献资料或者预实验，发现该载体在目标细胞模型中表达效率过低，或者会受到实验研究因素干扰时，再考虑其他表达活性更强的载体（如 pGL4.75），或自己构建所需的特殊启动子载体。
- 如果萤光素酶表达量及信号强度满足要求的前提下，可选择快速应答报告基因载体，即萤光素酶基因与 PEST 失稳序列的融合（图 A），该融合缩短了萤光素酶的细胞半衰期 ($t_{1/2}$)，防止了报告基因蛋白的积累，从而能够更快更灵敏的检测诱导反应（图 B & C），更容易区别阳性和阴性信号。



双萤光素酶报告基因载体列表

RE 克隆用骨架载体

骨架元件	报告基因缩写	稳转筛选 Marker	说明	产品	目录号
NanoLuc® 萤光素酶报告基因	Nluc	—	无快速应答功能, 可克隆已知或预测的启动子区域	pNL1.1 [Nluc] Vector	N1001
	NlucP	—	中度快速应答, 可克隆已知或预测的启动子区域, 报告基因在细胞内表达低于 Nluc	pNL1.2 [NlucP] Vector	N1011
	secNluc	—	分泌型 NanoLuc® 报告基因载体, 可克隆入目标启动子	pNL1.3 [secNluc] Vector	N1021
	Nluc	Hygro	潮霉素筛选 Marker, 可筛选稳转细胞系	pNL2.1 [Nluc/ Hygro] Vector	N1061
	NlucP	Hygro	潮霉素筛选 Marker, 中度快速应答, 可筛选稳转细胞系	pNL2.2 [NlucP/ Hygro] Vector	N1071
	secNluc	Hygro	分泌型萤光素酶, 潮霉素筛选 Marker, 可筛选稳转细胞系	pNL2.3 [secNluc/ Hygro] Vector	N1081
萤火虫萤光素酶报告基因	Luc2	—	无快速应答功能, 可克隆入目标启动子报告基因	pGL4.10[luc2] Vector	E6651
	Luc2P	—	中度快速应答, 可克隆入研究用启动子报告基因, 报告基因在细胞内表达低于 luc2, 高于 luc2CP	pGL4.11[luc2P] Vector	E6661
	Luc2CP	—	高度快速应答, 可克隆入研究用启动子报告基因, 报告基因在细胞内表达低于 luc2, luc2P	pGL4.12[luc2CP] Vector	E6671
	Luc2	Neo	新霉素筛选 Marker, 可筛选稳转细胞系	pGL4.17[luc2/ Neo] Vector	E6721
	Luc2	Hygro	潮霉素筛选 Marker, 可筛选稳转细胞系	pGL4.14[luc2/ Hygro] Vector	E6691
	Luc2	Puro	嘌呤霉素筛选 Marker, 可筛选稳转细胞系	pGL4.20[luc2/ Puro] Vector	E6751
海肾萤光素酶报告基因	hRluc	—	无快速应答功能, 报告基因在细胞内表达高	pGL4.70[hRluc] Vector	E6881

含 minP 或组成型启动子载体

报告基因	报告基因缩写	启动子缩写	说明	产品	目录号
NanoLuc® 萤光素酶报告基因	Nluc	minP	TATA box	pNL3.1 [Nluc/ minP] Vector	N1031
	NlucP	minP	TATA box	pNL3.2 [NlucP/ minP] Vector	N1041
	secNluc	minP	TATA box	pNL3.3 [secNluc/ minP] Vector	N1051
	Nluc	PGK	哺乳动物来源启动子, 强度适中	pNL1.1.PGK [Nluc/ PGK] Vector	N1441
	Nluc	TK	病毒来源组成型启动子, 强度低于 SV40 和 CMV	pNL1.1.TK[Nluc/ TK] Vector	N1501
	Nluc	CMV	病毒来源组成型启动子, 强度高于 SV40, TK, 在初次实验时的首选	pNL1.1.CMV[Nluc/CMV] Vector	N1091
	NlucP	CMV	病毒来源组成型启动子, 强度高于 SV40, TK, 中度快速应答	pNL3.2.CMV[NlucP/CMV] Vector	N1411
	secNluc	CMV	病毒来源组成型启动子, 强度高于 SV40, TK, 分泌型	pNL1.3.CMV[secNluc/ CMV] Vector	N1101
萤火虫萤光素酶报告基因	Luc2	PGK	哺乳动物来源启动子, 强度适中	pGL4.53[luc2/ PGK] Vector	E5011
	Luc2	minP	TATA box	pGL4.23 [luc2/ minP] Vector	E8411
	Luc2	TK	病毒来源组成型启动子, 强度低于 SV40 和 CMV	pGL4.54[luc2/ TK] Vector	E5061
	Luc2	SV40	病毒来源组成型启动子, 强度低于 CMV, 高于 TK	pGL4.13[luc2/ SV40] Vector	E6681
	Luc2	CMV	病毒来源组成型启动子, 强度高于 SV40, TK	pGL4.50[luc2/CMV/ Hygro] Vector	E1310
海肾萤光素酶报告基因	hRluc	TK	带有 TK 启动子的海肾萤光素酶报告基因载体, 在初次实验时的首选	pGL4.74[hRluc/ TK] Vector	E6921
	hRluc	SV40	带有 SV40 启动子的海肾萤光素酶报告基因载体, 强度低于 CMV, 高于 TK	pGL4.73[hRluc/ SV40] Vector	E6911
	hRluc	CMV	带有 CMV 启动子的海肾萤光素酶报告基因载体, 强度高于 TK 和 SV40	pGL4.75[hRluc/ CMV] Vector	E6931

预构建载体列表

商品化预构建信号通路元件载体，客户可根据实验需要直接选择相应载体购买，无需自行构建。

进行瞬时转染操作时，一般需设计阴性和阳性对照组，例如：需选择一个 *Renilla* Luciferase 的内参载体，还需选择 Firefly Luciferase 的空骨架载体（阴性对照组）及一个带有已知会受待测药物或其他物质影响的调控序列的 Firefly Luciferase 载体（阳性对照组）。

Response Element/Promoter 应答元件 / 启动子	Pathway 信号通路	Reporter Gene 报告基因	Vector Name 载体名称	目录号
Human PSA Promoter	Androgen Activation in Prostate Cancer	Firefly	pGL4[luc2P/PSA-long/Hygro] Vector	咨询 Promega
Cyclic AMP Response Element (CRE)	cAMP/PKA	NanoLuc [®]	pNL[NlucP/CRE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.29[luc2P/CRE/Hygro] Vector	E8471
NFAT Response Element	Calcium/Calcineurin	NanoLuc [®]	pNL[NlucP/NFAT-RE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.30[luc2P/NFAT-RE/Hygro] Vector	E8481
Human 2B6	Cytochrome 450/Drug Metabolism	Firefly	pGL4[luc2P/2B6/Hygro] Vector	咨询 Promega
Human 3A4		Firefly	pGL4[luc2P/3A4/Hygro] Vector	咨询 Promega
p53 Response Element (p53 RE)	DNA Damage/p53	NanoLuc [®]	pNL[NlucP/p53-RE/Hygro] Vector	咨询 Promega
		Firefly	pGL4[luc2P/p53-RE/Puro] Vector	咨询 Promega
		Firefly	pGL4.38[luc2P/p53 RE/Hygro] Vector	E3651
		<i>Renilla</i>	pGL4[hRlucP/p53-RE/Hygro] Vector	咨询 Promega
ATF4 Response Element (ATF4 ERSE)	Endoplasmic reticulum stress	NanoLuc [®]	pNL[NlucP/ATF4-RE/Hygro] Vector	咨询 Promega
		Firefly	pGL4[luc2P/ATF4-RE/Hygro] Vector	咨询 Promega
ATF6 Response Element (ATF6 ERSE)		NanoLuc [®]	pNL[NlucP/ATF6-RE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.39[luc2P/ATF6 RE/Hygro] Vector	E3661
Heat Shock Element (HSE)	Heat shock	Firefly	pGL4.41[luc2P/HSE/Hygro] Vector	E3751
Metal Regulatory Element (MRE)	Heavy metal stress	NanoLuc [®]	pNL[NlucP/MRE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.40[luc2P/MRE/Hygro] Vector	E4131
Gli Response Element (Gli)	Hedgehog	Firefly	pGL4[luc2P/Gli-RE/Hygro] Vector	咨询 Promega
Hypoxia Response Element (HRE)	Hypoxia	NanoLuc [®]	pNL[NlucP/HRE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.42[luc2P/HRE/Hygro] Vector	E4001
		<i>Renilla</i>	pGL4[hRlucP/HRE/Hygro] Vector	咨询 Promega
Human IL8 Promoter	IL1	Firefly	pGL4[luc2P/hIL8/Hygro] Vector	咨询 Promega
STAT5 Response Element	IL3	Firefly	pGL4.52[luc2P/STAT5 RE/Hygro]	E4651
sis-Inducible Element (SIE)	IL6	Firefly	pGL4.47[luc2P/SIE/Hygro] Vector	E4041
Human LCN2 (Lipocalin) Promoter	IL17	Firefly	pGL4[luc2P/LCN2/Hygro] Vector	咨询 Promega
Interferon Stimulated Response Element (ISRE)	JAK/STAT1/2 IFN α	NanoLuc [®]	pNL[NlucP/ISRE/Hygro] Vector	咨询 Promega
	INF- α	Firefly	pGL4.45[luc2P/ISRE/Hygro] Vector	E4141
IFN γ Activation Site (GAS) Response Element	JAK/STAT1 IFN γ	Firefly	pGL4[luc2P/GAS-RE/Hygro] Vector	咨询 Promega
		NanoLuc [®]	pNL[NlucP/GAS-RE/Hygro] Vector	咨询 Promega
Sis-Inducible Element (SIE) Response Element	AK/STAT3 IL6	NanoLuc [®]	pNL[NlucP/SIE/Hygro] Vector	咨询 Promega
IRF1 Response Element	JAK/STAT4 IL12	Firefly	pGL4[luc2P/STAT4-RE/Hygro] Vector	咨询 Promega
STAT5 Response Element	JAK/STAT5 IL3	NanoLuc [®]	pNL[NlucP/STAT5-RE/Hygro] Vector	咨询 Promega
Serum Response Element	MAP/ERK	Firefly	pGL4.33[luc2P/SRE/Hygro] Vector	E1340
AP1 Response Element (AP1 RE)	MAPK/JNK	NanoLuc [®]	pNL[NlucP/AP1-RE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.44[luc2P/AP1 RE/Hygro] Vector	E4111
		<i>Renilla</i>	pGL4[hRlucP/AP1-RE/Hygro] Vector	咨询 Promega
C/EBP Response Element (C/EBP)	Multiple	Firefly	pGL4[luc2P/C/EBP-RE/Hygro] Vector	咨询 Promega
		NanoLuc [®]	pNL[NlucP/C/EBP-RE/Hygro] Vector	咨询 Promega
Myc Response Element (Myc)	Myc	Firefly	pGL4[luc2P/Myc/Hygro] Vector	咨询 Promega
Nuclear Factor κ B Response Element	NF- κ B	NanoLuc [®]	pNL3.2.NF- κ B-RE[NlucP/NF- κ B-RE/Hygro] Vector	N1111
		Firefly	pGL4.32[luc2P/NF- κ B-RE/Hygro] Vector	E8491
NFAT Response Element	NFAT	Firefly	pGL4[luc2P/NFAT-RE/Puro] Vector	咨询 Promega
Human Early Growth Response Factor1 (EGR 1) Promoter	NGF	Firefly	pGL4[luc2P/hEGR1/Hygro] Vector	咨询 Promega
CBF1/RBPJk Response Element	Notch	Firefly	pGL4[luc2P/RBP-Jk-RE/Hygro] Vector	咨询 Promega

预构建载体列表

Response Element/Promoter 应答元件 / 启动子	Pathway 信号通路	Reporter Gene 报告基因	Vector Name 载体名称	目录号
Antioxidant Response Element (ARE)	Oxidative Stress	NanoLuc®	pNL[NlucP/ARE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.37[luc2P/ARE/Hygro] Vector	E3641
Myc/Max Response Element	PI3K/Akt/MAPK	NanoLuc®	pNL[NlucP/MycMax-RE/Hygro] Vector	咨询 Promega
Serum Response Element (SRE)	Ras/MEK1	NanoLuc®	pNL[NlucP/SRE/Hygro] Vector	咨询 Promega
Serum Response Factor (SRF) Response Element	RhoA (Gα12/13)	Firefly	pGL4[Luc2P/SRF-RE/Puro] Vector	咨询 Promega
		NanoLuc®	pNL[NlucP/SRF/Hygro] Vector	咨询 Promega
Serum Response Factor Response Element	RhoA	Firefly	pGL4.34[luc2P/SRF-RE/Hygro] Vector	E1350
Human G-CSF	STAT3	Firefly	pGL4[luc2P/GCSF/Hygro] Vector	咨询 Promega
Human IL2 Promoter	T Cell Activation IL1	Firefly	pGL4[luc2P/IL-2/Hygro] Vector	咨询 Promega
		Firefly	pGL4[luc2P/IL-2/Puro] Vector	咨询 Promega
SMAD3/SMAD4 Binding Element (SBE)	TGF-β	NanoLuc®	pNL[NlucP/SBE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.48[luc2P/SBE/Hygro] Vector	E3671
Human DNA-Binding Protein Inhibitor ID-1 Promoter	TGFβ/BMP	Firefly	pGL4[luc2P/hID1/Hygro] Vector	咨询 Promega
EGFR2 Promoter	VEGF2	Firefly	pGL4[Luc2P/hVEGFR2(-290 to +269)/Hygro] Vector	咨询 Promega
TCF-LEF Response Element (TCF-LEF RE)	Wnt	NanoLuc®	pNL[NlucP/TCF/LEF-RE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.49[luc2P/TCF-LEF RE/Hygro] Vector	E4611
Xenobiotic Responsive Element (XRE)	Xenobiotic Stress	NanoLuc®	pNL[NlucP/XRE/Hygro] Vector	咨询 Promega
		Firefly	pGL4.43[luc2P/XRE/Hygro] Vector	E4121

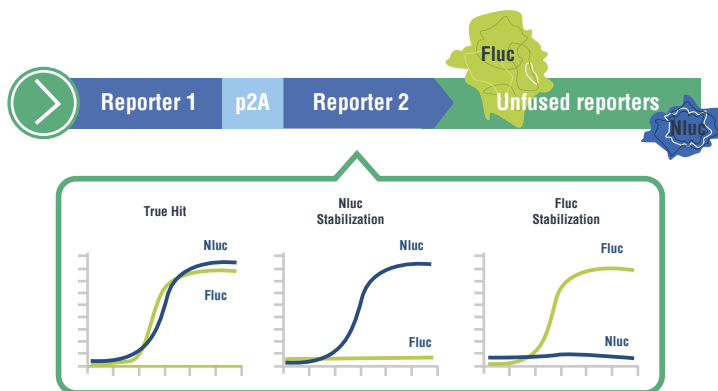
Coincidence 报告基因载体

两个报告基因，一个转录本

Coincidence 报告基因检测原理

通常，待测化合物与萤光素酶报告基因的相互作用往往也会导致一定的假阳性结果。此载体系统正是为此而设计。

原理：P2A 序列启动了核糖体跳跃，同一启动子可调控获得两种非融合酶的表达，且两种萤光素酶与化合物相互作用不同。因此，当用于高通量化合物筛选时，实现了与其中一种萤光素酶直接作用而导致的假阳性化合物能够与同时引起两种萤光素酶反应的真正有效化合物区别开来。



特点：

- pNLCol 载体能从相同的 mRNA 转录本中同时表达 Firefly 萤光素酶 (luc2) 和带有一个 PEST 去稳结构域的 NanoLuc® 萤光素酶 (NlucP)。通过来源于猪捷申病毒 -1 (porcine teschovirus-1) 的 P2A 序列能够分别检测两种萤光素酶的表达。
- pNLCol 载体设计是与 Nano-Glo® Dual-Luciferase® Reporter 检测系统一起使用，该系统以简单的“加样 - 读数 - 加样 - 读数”模式对同一样品中的 Firefly 和 NanoLuc® 萤光素酶活性进行顺序检测。

Promoter 启动子	Illustration 说明	Reporter Gene 报告基因	Vector Name 载体名称	目录号
--	骨架载体，包含多克隆位点区域，可插入感兴趣的启动子序列。	Firefly/NanoLuc®	pNLCol1[luc2-P2A-NlucP/Hygro] Vector	N1461
minP	包含用于在最小启动子 (minP) 上游插入启动子元件的多克隆位点区域。	Firefly/NanoLuc®	pNLCol2[luc2-P2A-NlucP/minP/Hygro] Vector	N1471
CMV	启动 Firefly/NanoLuc® 两个报告基因的高表达；可作为组成型表达的阳性对照载体。	Firefly/NanoLuc®	pNLCol3[luc2-P2A-NlucP/CMV/Hygro] Vector	N1481
PGK	启动 Firefly/NanoLuc® 两个报告基因的中度表达；可作为组成型表达的阳性对照载体。	Firefly/NanoLuc®	pNLCol4[luc2-P2A-NlucP/PGK/Hygro] Vector	N1491

双萤光素酶报告基因检测试剂

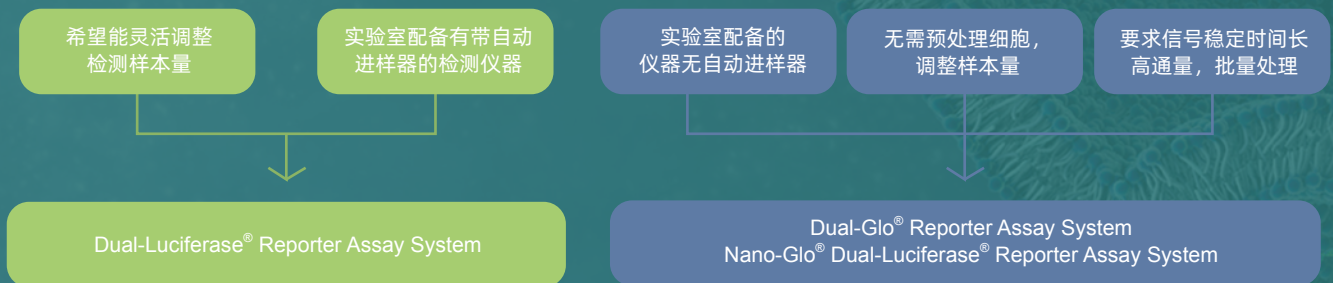
Promega 双萤光素酶报告基因专利检测系统可在同一个样品孔中先后检测 2 个不同的萤光素酶发光信号。试剂系统分为非均质法和均质法两种。客户可根据实验的需求和仪器条件进行选择。

• 如何选择试剂？

① 根据载体组合选择

载体组合 (主报告基因 / 内参报告基因)	产品	目录号	规格 (96 孔板)	特点
Firefly/Renilla	Dual-Luciferase [®] Reporter Assay System (缩写 DLR)	E1910	10ml	<ul style="list-style-type: none"> 非均质法，先去除上清，PBS 洗涤，裂解细胞，再加入底物试剂检测 改良型闪光法试剂，萤火虫萤光素酶光信号半衰期 10min 检测体积可灵活调整 如果检测样本较多，建议使用自动进样器加样
		E1960	10x10ml	
		E1980	100ml	
	Dual-Glo [®] Reporter Assay System	E2920	10ml	
		E2940	100ml	
		E2980	10X100ml	
NanoLuc [®] /Firefly 或 Firefly/NanoLuc [®]	Nano-Glo [®] Dual-Luciferase [®] Reporter Assay System	N1610	10ml	<ul style="list-style-type: none"> 均质法，无需去除上清和单独裂解细胞，两步加入试剂，操作更简便 光信号稳定时间更长，信号稳定时间长达 2h 排枪操作即可，无需自动进样器 但如果检测样本较多，仍建议采用自动进样器
		N1620	100ml	
		N1630	10X10ml	
		N1650	10X100ml	

② 均质法与非均质法的选择



• 举例比较均质法 (Dual-Glo[®]) 与非均质法 (Dual-Luciferase[®]) 操作流程：



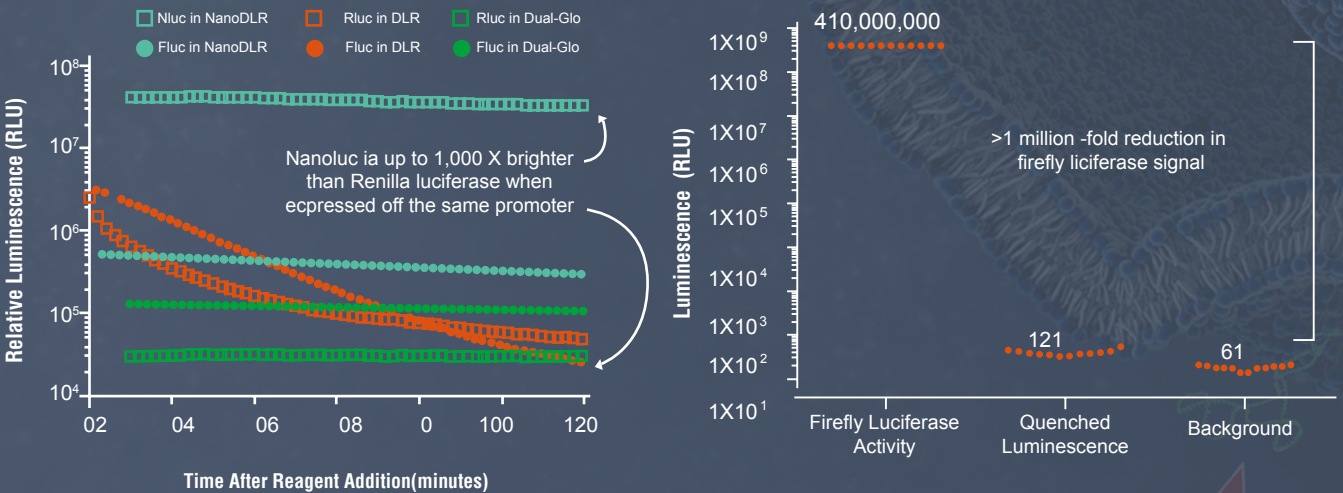
双报告基因检测全新升级版

• Nano-Glo® Dual-Luciferase® 双报告基因检测系统

Nano-Glo® Dual-Luciferase® Reporter (NanoDLR™) Assay System 是 Promega 全新推出的一款双报告基因检测系统，系统提供均质法检测试剂，能够在同一个样品中先后检测萤火虫萤光素酶 (*Photinus pyralis*) 和 NanoLuc® 萤光素酶 (Nluc) 的活性。先以 ONE-Glo™ EX Luciferase Assay Reagent 检测萤火虫萤光素酶活性，再以 NanoDLR™ Stop & Glo® Reagent 淬灭萤火虫萤光素酶信号，同时 NanoLuc® 的底物 Furimazine 与 NanoLuc® 反应产生光信号。这种“加样 - 检测 - 加样 - 检测”的模式能够产生稳定的辉光信号。

新系统亮点

- 双报告基因检测 - 萤火虫和 NanoLuc® 萤光素酶
- 辉光型检测信号 - 信号半衰期长，适合高通量检测
- 提供双报告基因检测所需的基础载体 - 主报告基因载体和内参载体
- 可在原有双报告基因系统（萤火虫 / 海肾）上灵活升级 - 仅需构建极少载体或使用商品化对照载体（含 NLuc）即可转换新系统
- 萤火虫和 NanoLuc® 萤光素酶任一均可作为主报告基因，也可将二者都作为主报告基因
- Nluc 信号更亮，所用载体量更少



如图所示，我们比较了 TK-Rluc (*Renilla*) : TK-Fluc (*Firefly*) : 载体 DNA 或者是 TK-Nluc (NanoLuc) : TK-Fluc (*Firefly*) : 载体 DNA 以 1:1:8 的比例转染 HEK293 细胞，通过 NanoDLR™, Dual-Glo® 或者 Dual-Luciferase® Reporter Assay System 测得的发光信号。

图中数据显示了在 NanoDLR™ 检测系统中萤火虫萤光素酶信号的有效淬灭。NanoDLR™ Stop & Glo® Reagent 可以将极高的 Fluc 信号降低至接近本底水平，这将十分有利于信号分离并减少变异性发生，且无需牺牲两个具有明亮发光信号酶的灵敏度。

相关产品

无内毒素质粒提取—PureYield™ 去内毒素质粒纯化系统!

- 系统包含独特的内毒素清除洗脱液，用于去除纯化质粒 DNA 中的蛋白质、RNA 和内毒素的污染。
- 使用新型硅基质膜，无需异丙醇沉淀或长时间的离心步骤，使操作更简单，纯化产量更高。

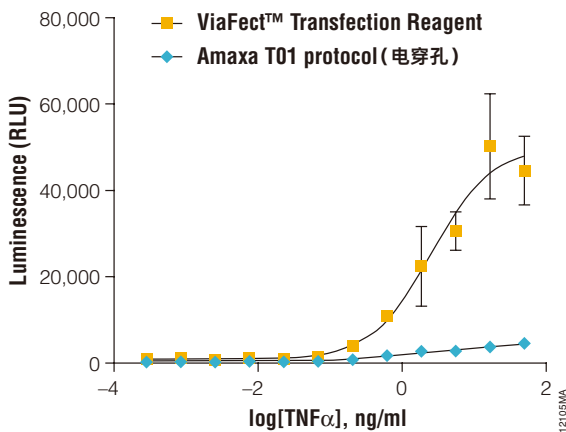
名称	产品特点	规格	目录号
PureYield™ Plasmid Miniprep System	<ul style="list-style-type: none"> • 仅需 10 分钟就能从 600µl 到 3ml 细菌培养物中纯化出高达 15µg 的质粒 DNA • 无需异丙醇沉淀或长时间的离心步骤 • 可使用离心或真空法操作 	100 次	A1223
		250 次	A1222
PureYield™ Plasmid Maxiprep System	<ul style="list-style-type: none"> • 从 250ml 培养液中提取高达 1mg 转染级别质粒 • 基于真空的方案 • 高纯度浓缩质粒 DNA • 没有冗长的重力滴灌注或洗脱后醇沉淀步骤 	10 preps	A2392
		25 preps	A2393

高效低毒转染 -ViaFect™ 转染试剂

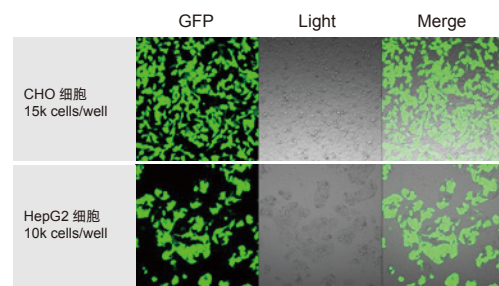
下图：以 ViaFect™ 转染试剂建立造血细胞的细胞因子信号通路模型。以 ViaFect™ 转染试剂或电穿孔将 pGL4.32[*luc2P*/NF-κB-RE/Hygro] Vector (一种带有 NF-κB 应答元件的萤光素酶报告基因载体) 瞬时转染入 TF-1 cells。然后以 TNFα 刺激细胞 6 小时，再以 Bio-Glo™ Luciferase Assay System (用于 Biologics 检测) 检测萤光素酶表达。



在难转染的悬浮细胞中建立细胞监测模型



转染效果举例



*注：反向转染指先在培养板孔中加入 DNA：转染试剂混合物，再加入细胞进行转染的方式。

产品功能	产品	规格	目录号	说明
将载体转染入细胞	ViaFect™ Transfection Reagent	0.2ml	E4983	<ul style="list-style-type: none"> • 创新型配方转染试剂； • 性价比更高； • 超低细胞毒性； • 对于难转染的多种细胞效果更佳； • 操作简单，无需更换培养基。
		0.75ml	E4981	
		2x 0.75ml	E4982	

GloMax[®] Discover 多功能检测仪

高效，易用的多功能检测仪：用于化学发光，荧光，吸收光，BRET 和 FRET

- **使用简单**

选择预置的 Promega 程序或编辑您自己需要的程序。数据可输出至网络，云，LIMS 或其他服务器。容易实现化学发光，荧光，UV-可见吸收光，BRET 和 FRET 检测。

- **可实现自动化**

可与您的自动化工作流程整合实现高通量检测，或者整合进您的 LIMS 数据管理系统。

- **超高灵敏度**

动态范围宽，灵敏度高，孔间交叉干扰小，使得您的数据更加可信。

- **手动干扰最小化**

自动转换滤片，轻松实现多重检测。

- **节省人力**

预置的程序可节省优化程序的时间。

- **直观的软件界面**

触摸屏可实现编辑程序，整合自动化平台，输出数据。



监测仪器	目录号	包装量
GloMax [®] Discover System	GM3000	1 台

萤光素酶报告基因技术应用文献

应用	文献
病毒研究	<p>Chao Zhang, Yifan Wang, Yuanfei Zhu, et al. (2021) Development and structural basis of a two-Mab cocktail for treating SARS-CoV-2 infections. <i>Nat Commun.</i> 12(1):264 [IF: 14.919]</p> <p>Chao Zhang, Cong Xu, Wenlong Dai, et al. (2021) Functional and structural characterization of a two-MAb cocktail for delayed treatment of enterovirus D68 infections. <i>Nat Commun.</i> 12(1):2904 [IF: 14.919]</p> <p>Zhongcheng Zhou, Peng Du, Meixing Yu, et al. (2021) Assessment of infectivity and the impact on neutralizing activity of immune sera of the COVID-19 variant, CAL.20C. <i>Signal Transduction and Targeted Therapy.</i> 6:285 [IF: 18.187]</p>
疾病研究 (糖尿病肾病, 皮肤瘤, 黑素瘤等)	<p>Huiming Liu, Rui Yan, Luqun Liang, et al. (2021) The role of CDX2 in renal tubular lesions during diabetic kidney disease. <i>Aging (Albany NY).</i> 13(5): 6782-6803 [IF: 5.682]</p> <p>Hui-Na Wang, Kunmei Ji, Li-Na Zhang, et al. (2021) Inhibition of c-Fos expression attenuates IgE-mediated mast cell activation and allergic inflammation by counteracting an inhibitory AP1/Egr1/IL-4 axis. <i>Journal of Translational Medicine.</i> 6:285 [IF: 5.531]</p> <p>Yeye Guo, Xu Zhang, Weiqi Zeng, et al. (2020) TRAF6 Activates Fibroblasts to Cancer-Associated Fibroblasts through FGF19 in Tumor Microenvironment to Benefit the Malignant Phenotype of Melanoma Cells. <i>J Invest Dermatol.</i> 140(11):2268-2279 [IF: 8.551]</p> <p>Xu Zhang, Lisha Wu, Ta Xiao, et al. (2018) TRAF6 regulates EGF-induced cell transformation and cSCC malignant phenotype through CD147/EGFR. <i>Oncogenesis.</i> 7(2): 17 [IF: 7.485]</p>
代谢研究	<p>Chi Qianru, Hu Xueyuan, Zhao Bing, et al. (2021) Regulation of H2S-induced necroptosis and inflammation in broiler bursa of Fabricius by the miR-15b-5p/TGFBR3 axis and the involvement of oxidative stress in this process. <i>Journal of Hazardous Materials.</i> 406:124682 [IF: 10.588]</p>
调控元件功能研究 (启动子, 增强子等)	<p>Andrienas K.K., Ramlall V., Kurland J., et al. (2018). DNA-binding landscape of IRF3, IRF5 and IRF7 dimers: implications for dimer-specific gene regulation. <i>Nucleic Acids Res.</i> 46(5):2509-2520 [IF: 16.971]</p> <p>Cacace A, Sboarina M, Vazeille T, et al. (2017). Glutamine activates STAT3 to control cancer cell proliferation independently of glutamine metabolism. <i>Oncogene.</i> 36(15):2074-2084 [IF: 9.867]</p>
分子马达	<p>Xia Xu, Jinbo Fei, Youqian Xu et al. (2021) Boric Acid-Fueled ATP Synthesis by FoF1 ATP Synthase Reconstituted in a Supramolecular Architecture. <i>Angewandte Chemie International Edition.</i> 60(14):7617-7620 [IF: 15.336]</p>

萤光素酶报告基因技术应用文献 (接上页)

应用	文献
转录因子活性或细胞信号通路传导	<p>Kankia, I.H., Khalil, H.S., Langdon, S.P., et al. (2017). NRF2 Regulates HER1 Signaling Pathway to Modulate the Sensitivity of Ovarian Cancer Cells to Lapatinib and Erlotinib. <i>Oxid. Med. Cell Longev.</i> 2017:1864578 [IF: 6.543]</p> <p>Lynch RA, Etschin J, Battle TE, et al. (2007). A small-molecule enhancer of signal transducer and activator of transcription 1 transcriptional activity accentuates the antiproliferative effects of IFN-gamma in human cancer cells. <i>Cancer Res.</i> 67(3):1254-61 [IF: 12.701]</p>
膜受体研究	<p>Kimble AJ, Soundararajan M, Hutsell SQ, et al. (2009). Structural Determinants of G-protein Subunit Selectivity by Regulator of G-protein Signaling 2 (RGS2). <i>J. Biol. Chem.</i> 284(29):19402-11 [IF: 5.157]</p>
双杂交法蛋白相互作用	<p>Traustadóttir GÁ, Jensen CH, Garcia Ramirez JJ, et al. (2017). The non-canonical NOTCH1 ligand Delta-like 1 homolog (DLK1) self interacts in mammals. <i>Int J Biol Macromol.</i> 97:460-467 [IF: 6.953]</p>
siRNA/miRNA	<p>Hatice S Kaya-Okur, Steven J Wu, Christine A Codomo, et al. (2019) CUT&Tag for efficient epigenomic profiling of small samples and single cells. <i>Nat Commun.</i> 10(1):1930 [IF: 14.919]</p> <p>Soldà G, Robusto M, Primignani P, et al. (2012). A novel mutation within the MIR96 gene causes non-syndromic inherited hearing loss in an Italian family by altering pre-miRNA processing. <i>Hum Mol Genet.</i> 21(3):577-85 [IF: 6.15]</p>
植物研究	<p>Wan Zhao, Jinting Yu, Feng Jiang, et al. (2021) Coordination between terminal variation of the viral genome and insect microRNAs regulates rice stripe virus replication in insect vectors. <i>PLoS Pathog.</i> 17(3):e1009424 [IF: 6.823]</p> <p>Hong Liu, Huifang, Chenyang Hao, et al. (2020) TaDA1, a conserved negative regulator of kernel size, has an additive effect with TaGW2 in common wheat (<i>Triticum aestivum</i> L.) <i>Plant Biotechnol J.</i> 18(5):1330-1342 [IF: 9.803]</p> <p>Rahmani, F, Hummel, M, Schuurmans, J, et al. (2009). Sucrose control of translation mediated by an upstream open reading frame-encoded peptide. <i>Plant Physiol.</i> 150(3):1356-67 [IF: 8.34]</p> <p>Dunoyer P, Lecellier CH, Parizotto EA, et al. (2004). Probing the microRNA and small interfering RNA pathways with virus-encoded suppressors of RNA silencing. <i>Plant Cell.</i> 16(5):1235-50 [IF: 11.277]</p>

萤光素酶报告基因载体

载体	报告基因	启动子	PEST 序列	筛选标记	目录号
NanoLuc® 报告基因载体					
无启动子载体					
pNL1.1 [Nluc] Vector	Nluc	-	-	-	N1001
pNL1.2 [NlucP] Vector	NlucP	-	hPEST	-	N1011
pNL1.3 [secNluc] Vector	secNluc	-	-	-	N1021
pNL2.1 [Nluc/Hygro] Vector	Nluc	-	-	Hygro	N1061
pNL2.2 [NlucP/Hygro] Vector	NlucP	-	hPEST	Hygro	N1071
pNL2.3 [secNluc/Hygro] Vector	secNluc	-	-	Hygro	N1081
含 minP 载体					
pNL3.1 [Nluc/minP] Vector	Nluc	minP	-	-	N1031
pNL3.2 [NlucP/minP] Vector	NlucP	minP	hPEST	-	N1041
pNL3.3 [secNluc/minP] Vector	secNluc	minP	-	-	N1051
pNL[NlucP/minP/Hygro] Vector	NlucP	minP	hPEST	Hygro	咨询 Promega
内参报告基因载体					
pNL1.1.CMV [Nluc/CMV] Vector	Nluc	CMV	-	-	N1091
pNL1.1.PGK[Nluc/PGK] Vector	Nluc	PGK	-	-	N1441
pNL1.1.TK[Nluc/TK]Vector	Nluc	TK	-	-	N1501
pNL1.3.CMV [secNluc/CMV] Vector	secNluc	CMV	-	-	N1101
pNL3.2.CMV[NlucP/CMV] Vector	NlucP	CMV	hPEST	-	N1411
Coincidence reporter 载体					
pNLCol1[luc2-P2A-NlucP/Hygro] Vector	Luc2, NlucP	-	hPEST	Hygro	N1461
pNLCol2[luc2-P2A-NlucP/minP/Hygro] Vector	Luc2, NlucP	-	hPEST	Hygro	N1471
pNLCol3[luc2-P2A-NlucP/CMV/Hygro] Vector	Luc2, NlucP	-	hPEST	Hygro	N1481
pNLCol4[luc2-P2A-NlucP/PGK/Hygro] Vector	Luc2, NlucP	-	hPEST	Hygro	N1491
Firefly 报告基因载体					
无启动子载体					
pGL4.10[luc2] Vector	Luc2	-	-	-	E6651
pGL4.11[luc2P] Vector	Luc2P	-	hPEST	-	E6661
pGL4.12[luc2CP] Vector	Luc2CP	-	hPEST	-	E6671

萤光素酶报告基因载体 (接上页)

载体	报告基因	启动子	PEST 序列	筛选标记	目录号
Firefly 报告基因载体					
无启动子载体					
pGL4.14[luc2/ Hygro] Vector	Luc2	-	-	Hygro	E6691
pGL4.15[luc2P/ Hygro] Vector	Luc2P	-	hPEST	Hygro	E6701
pGL4.16[luc2CP/ Hygro] Vector	Luc2CP	-	hPEST	Hygro	E6711
pGL4.17[luc2/ Neo] Vector	Luc2	-	-	Neo	E6721
pGL4.18[luc2P/ Neo] Vector	Luc2P	-	hPEST	Neo	E6731
pGL4.19[luc2CP/ Neo] Vector	Luc2CP	-	hPEST	Neo	E6741
pGL4.20[luc2/ Puro] Vector	Luc2	-	-	Puro	E6751
pGL4.21[luc2P/ Puro] Vector	Luc2P	-	hPEST	Puro	E6761
pGL4.22[luc2CP/ Puro] Vector	Luc2CP	-	hPEST	Puro	E6771
含 minP 载体					
pGL4.23[luc2/minP]Vector	Luc2	minP	-	-	E8411
pGL4.24[luc2P/minP]Vector	Luc2P	minP	hPEST	-	E8421
pGL4.25[luc2CP/minP]Vector	Luc2CP	minP	hPEST	-	E8431
pGL4.26[luc2/minP/Hygro]Vector	Luc2	minP	-	Hygro	E8441
pGL4.27[luc2P/minP/Hygro]Vector	Luc2P	minP	hPEST	Hygro	E8451
pGL4.28[luc2CP/minP/Hygro]Vector	Luc2CP	minP	hPEST	Hygro	E8461
内参报告基因载体					
pGL4.13[luc2/ SV40] Vector	Luc2	SV40	-	-	E6681
pGL4.50[luc2/CMV/ Hygro] Vector	Luc2	CMV	-	Hygro	E1310
pGL4.51[luc2/CMV/ Neo] Vector	Luc2	CMV	-	Neo	E1320
pGL4.53[luc2/ PGK] Vector	Luc2	PGK	-	-	E5011
pGL4.54[luc2/ TK] Vector	Luc2	TK	-	-	E5061
融合载体					
pSP-luc+NF Fusion Vector (N-Term)	Luc+	SP6 (sense); T7 (antisense)	-	-	E4471

萤光素酶报告基因载体 (接上页)

载体	报告基因	启动子	PEST 序列	筛选标记	目录号
Renilla 报告基因载体					
无启动子载体					
pGL4.70[hRluc] Vector	hRluc	-	-	-	E6881
pGL4.82[hRluc/Puro] Vector	hRluc	-	-	Puro	E7501
pGL4.83[hRlucP/Puro] Vector	hRlucP	-	hPEST	Puro	E7511
pGL4.84[hRlucCP/Puro] Vector	hRlucCP	-	hPEST	Puro	E7521
pRL-null Vector	Rluc**	-	-	-	E2271
含 minP 载体					
pGL4[hRlucP/minP/Hygro]	hRlucP	minP	hPEST	Hygro	咨询 Promega
内参报告基因载体					
pGL4.73[hRluc/ SV40] Vector	hRluc	SV40	-	-	E6911
pGL4.74[hRluc/ TK] Vector	hRluc	TK	-	-	E6921
pGL4.75[hRluc/ CMV] Vector	hRluc	CMV	-	-	E6931
pRL-CMV Vector	Rluc**	CMV	-	-	E2261
pRL-SV40 Vector	Rluc**	SV40	-	-	E2231
pRL-TK Vector	Rluc**	TK	-	-	E2241
RNA 干扰					
pmirGLO Vector	luc2, hRluc	PGK, SV40	-	Neo	E1330
pmirNanoGlo Vector	luc2, NlucP	SV40, PGK	PEST	Blasticidin	咨询 Promega
psiCHECK™-2 Vector	hRluc, hluc+	SV40, HSV-TK	-	-	C8021

** wild-type *Renilla* luciferase

萤光素酶报告基因载体 (接上页)

Pathway 信号通路	Transcription Factor 转录因子	Response Element/Promoter 应答元件 / 启动子	Reporter Gene 报告基因	Vector Name 载体名称	Catlog 目录号
GPCR 信号转导与细胞应激反应通路载体					
Oxidative Stress	Nrf2	Antioxidant Response Element (ARE)	NanoLuc [®]	pNL[NlucP/ARE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.37[luc2P/ARE/Hygro] Vector	E3641
cAMP/PKA	CREB	Cyclic AMP Response Element (CRE)	NanoLuc [®]	pNL[NlucP/CRE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.29[luc2P/CRE/Hygro] Vector	E8471
Calcium/Calcineurin	NFAT	NFAT Response Element	NanoLuc [®]	pNL[NlucP/NFAT-RE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.30[luc2P/NFAT-RE/Hygro] Vector	E8481
NF-κB	NFκB	Nuclear Factor κB Response Element	NanoLuc [®]	pNL3.2.NF-κB-RE[NlucP/NF-κB-RE/Hygro] Vector	N1111
			Firefly	pGL4.32[luc2P/NF-κB-RE/Hygro] Vector	E8491
RhoA (Gα12/13)	SRF	Serum Response Factor (SRF) Response Element	NanoLuc [®]	pNL[NlucP/SRF/Hygro] Vector	咨询 Promega
			Firefly	pGL4[luc2P/SRF-RE/Puro] Vector	咨询 Promega
RhoA	SRF	Serum Response Factor Response Element	Firefly	pGL4.34[luc2P/SRF-RE/Hygro] Vector	E1350
DNA Damage/p53	p53	p53 Response Element (p53 RE)	NanoLuc [®]	pNL[NLucP/p53-RE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.38[luc2P/p53 RE/Hygro] Vector	E3651
			Firefly	pGL4[luc2P/p53-RE/Puro] Vector	咨询 Promega
			<i>Renilla</i>	pGL4[hRlucP/p53-RE/Hygro] Vector	咨询 Promega
Endoplasmic reticulum stress	ATF4	ATF4 Response Element (ATF4 ERSE)	NanoLuc [®]	pNL[NlucP/ATF4-RE/Hygro] Vector	咨询 Promega
			Firefly	pGL4[luc2P/ATF4-RE/Hygro] Vector	咨询 Promega
Endoplasmic reticulum stress	ATF6	ATF6 Response Element (ATF6 ERSE)	NanoLuc [®]	pNL[NlucP/ATF6-RE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.39[luc2P/ATF6 RE/Hygro] Vector	E3661
Heat shock	HSF1	Heat Shock Element (HSE)	Firefly	pGL4.41[luc2P/HSE/Hygro] Vector	E3751
Hypoxia	HIF1 α	Hypoxia Response Element (HRE)	NanoLuc [®]	pNL[NlucP/HRE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.42[luc2P/HRE/Hygro] Vector	E4001
			<i>Renilla</i>	pGL4[hRlucP/HRE/Hygro] Vector	咨询 Promega
Xenobiotic Stress	AhR	Xenobiotic Responsive Element (XRE)	NanoLuc [®]	pNL[NlucP/XRE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.43[luc2P/XRE/Hygro] Vector	E4121
MAPK/JNK	AP1	AP1 Response Element (AP1 RE)	NanoLuc [®]	pNL[NlucP/AP1-RE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.44[luc2P/AP1 RE/Hygro] Vector	E4111
			<i>Renilla</i>	pGL4[hRlucP/AP1-RE/Hygro] Vector	咨询 Promega
PI3K/Akt/MAPK	Myc/Max	Myc/Max Response Element	NanoLuc [®]	pNL[NlucP/MycMax-RE/Hygro] Vector	咨询 Promega
Myc		Myc Response Element (Myc)	Firefly	pGL4[luc2P/Myc/Hygro] Vector	咨询 Promega
Androgen Activation in Prostate Cancer	-	Human PSA Promoter	Firefly	pGL4[luc2P/PSA-long/Hygro] Vector	咨询 Promega
Cytochrome 450/Drug Metabolism	-	Human 2B6	Firefly	pGL4[luc2P/2B6/Hygro] Vector	咨询 Promega
	-	Human 3A4	Firefly	pGL4[luc2P/3A4/Hygro] Vector	咨询 Promega
MAP/ERK	Elk-1/SRF	Serum Response Element (SRE)	Firefly	pGL4.33[luc2P/SRE/Hygro] Vector	E1340
Heavy metal stress	MTF1	Metal Regulatory Element (MRE)	NanoLuc [®]	pNL[NlucP/MRE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.40[luc2P/MRE/Hygro] Vector	E4131
Ras/MEK1	-	Serum Response Element (SRE)	NanoLuc [®]	pNL[NlucP/SRE/Hygro] Vector	咨询 Promega

萤光素酶报告基因载体 (接上页)

Pathway 信号通路	Transcription Factor 转录因子	Response Element/Promoter 应答元件 / 启动子	Reporter Gene 报告基因	Vector Name 载体名称	Catlog 目录号
细胞因子信号通路载体					
JAK/STAT1/2 IFN α	STAT1: STAT2	Interferon Stimulated Response Element (ISRE)	NanoLuc [®]	pNL[NlucP/ISRE/Hygro] Vector	咨询 Promega
INF- α			Firefly	pGL4.45[luc2P/ISRE/Hygro] Vector	E4141
JAK/STAT3 IL6	STAT3: STAT3	Sis-Inducible Element (SIE) Response Element	NanoLuc [®]	pNL[NlucP/SIE/Hygro] Vector	咨询 Promega
IL6			Firefly	pGL4.47[luc2P/SIE/Hygro] Vector	E4041
TGF- β	SMAD3 : SMAD4	SMAD3/SMAD4 Binding Element (SBE)	NanoLuc [®]	pNL[NlucP/SBE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.48[luc2P/SBE/Hygro] Vector	E3671
TGF β /BMP	-	Human DNA-Binding Protein Inhibitor ID-1 Promoter	Firefly	pGL4[luc2P/hID1/Hygro] Vector	咨询 Promega
Wnt	TCF-LEF	TCF-LEF Response Element (TCF-LEF RE)	NanoLuc [®]	pNL[NlucP/TCF/LEF-RE/Hygro] Vector	咨询 Promega
			Firefly	pGL4.49[luc2P/TCF-LEF RE/Hygro] Vector	E4611
JAK/STAT5 IL3	STAT5: STAT5	STAT5 Response Element	NanoLuc [®]	pNL[NlucP/STAT5-RE/Hygro] Vector	咨询 Promega
IL3			Firefly	pGL4.52[luc2P/STAT5 RE/Hygro]	E4651
IL1	-	Human IL8 Promoter	Firefly	pGL4[luc2P/hIL8/Hygro] Vector	咨询 Promega
T Cell Activation IL1	-	Human IL2 Promoter	Firefly	pGL4[luc2P/IL-2/Hygro] Vector	咨询 Promega
			Firefly	pGL4[luc2P/IL-2/Puro] Vector	咨询 Promega
IL17	-	Human LCN2 (Lipocalin) Promoter	Firefly	pGL4[luc2P/LCN2/Hygro] Vector	咨询 Promega
JAK/STAT1 IFN γ	IFN γ Activation Site	IFN γ Activation Site (GAS) Response Element	NanoLuc [®]	pNL[NlucP/GAS-RE/Hygro] Vector	咨询 Promega
			Firefly	pGL4[luc2P/GAS-RE/Hygro] Vector	咨询 Promega
JAK/STAT4 IL12	Response Element	IRF1 Response Element	Firefly	pGL4[luc2P/STAT4-RE/Hygro] Vector	咨询 Promega
Hedgehog	Gli	Gli Response Element (Gli-RE)	Firefly	pGL4[luc2P/Gli-RE/Hygro] Vector	咨询 Promega
Multiple	C/EBP	C/EBP Response Element (C/EBP)	NanoLuc [®]	pNL[NlucP/C/EBP-RE/Hygro] Vector	咨询 Promega
			Firefly	pGL4[luc2P/C/EBP-RE/Hygro] Vector	咨询 Promega
STAT3	-	Human G-CSF	Firefly	pGL4[luc2P/GCSF/Hygro] Vector	咨询 Promega
VEGF2	-	hVEGFR2 Promoter	Firefly	pGL4[Luc2P/hVEGFR2(-290 to +269)/Hygro] Vector	咨询 Promega
NGF	-	Human Early Growth Response Factor1 (EGR 1) Promoter	Firefly	pGL4[luc2P/hEGR1/Hygro] Vector	咨询 Promega
Notch	-	CBF1/RBPJk Response Element	Firefly	pGL4[luc2P/RBP-Jk-RE/Hygro] Vector	咨询 Promega
NFAT	-	NFAT Response Element	Firefly	pGL4[luc2P/NFAT-RE/Puro] Vector	咨询 Promega

萤光素酶报告基因检测

产品	规格	目录号
NanoLuc® 报告基因检测		
Nano-Glo® Luciferase Assay	10 ml	N1110
Nano-Glo® Luciferase Assay	100 ml	N1120
Nano-Glo® Luciferase Assay	10 x 10 ml	N1130
Nano-Glo® Luciferase Assay	10 x 100 ml	N1150
Firefly 报告基因检测		
Luciferase Assay Reagent	100 ml	E1483
Luciferase Assay System	100 assays	E1500
Luciferase Assay System, 10-Pack	1000 assays	E1501
Luciferase Assay System with Reporter Lysis Buffer	100 assays	E4030
Luciferase Assay System Freezer Pack	1000 assays	E4530
Luciferase 1000 Assay System	1000 assays	E4550
Luciferase Cell Culture Lysis 5X Reagent	30 ml	E1531
Reporter Lysis 5X Buffer	30 ml	E3971
Bright-Glo™ Luciferase Assay System	10 ml	E2610
Bright-Glo™ Luciferase Assay System	100 ml	E2620
Bright-Glo™ Luciferase Assay System	10 x 100 ml	E2650
ONE Glo™ Luciferase Assay System	10 ml	E6110
ONE Glo™ Luciferase Assay System	100 ml	E6120
ONE Glo™ Luciferase Assay System	1 L	E6130
ONE-Glo™ EX Luciferase Assay System	10 ml	E8110
ONE-Glo™ EX Luciferase Assay System	100 ml	E8120
ONE-Glo™ EX Luciferase Assay System	10 x 10 ml	E8130
ONE-Glo™ EX Luciferase Assay System	10 x 100 ml	E8150
Steady-Glo® Luciferase Assay System	10 ml	E2510
Steady-Glo® Luciferase Assay System	100 ml	E2520
Steady-Glo® Luciferase Assay System	10 x 100 ml	E2550
Renilla 报告基因检测		
Renilla Luciferase Assay System	100 assays	E2810
Renilla Luciferase Assay System	1000 assays	E2820
Renilla-Glo® Luciferase Assay System	10 ml	E2710
Renilla-Glo® Luciferase Assay System	100 ml	E2720
Renilla-Glo® Luciferase Assay System	10 x 100 ml	E2750

萤光素酶报告基因检测 (接上页)

产品	规格	目录号
NanoLuc® + Firefly 萤光素酶组成的双报告基因检测		
Nano-Glo® Dual-Luciferase® Reporter Assay System	10 ml	N1610
Nano-Glo® Dual-Luciferase® Reporter Assay System	100 ml	N1620
Nano-Glo® Dual-Luciferase® Reporter Assay System	10 x 10 ml	N1630
Nano-Glo® Dual-Luciferase® Reporter Assay System	10 x 100 ml	N1650
Nano-Glo® Dual-Luciferase® Reporter Assay/pNL1.1.TK Bundle	1 each	N1521
Nano-Glo® Dual-Luciferase® Reporter Assay/pNL1.1.PGK Bundle	1 each	N1531
Nano-Glo® Dual-Luciferase® Reporter Assay/pGL4.54[luc2/TK] Bundle	1 each	N1541
Nano-Glo® Dual-Luciferase® Reporter Assay/pGL4.53[luc2/PGK]Bundle	1 each	N1551
Passive Lysis 5X Buffer	30ml	E1941
Firefly + Renilla 萤光素酶组成的双报告基因检测		
Dual-Luciferase® Reporter Assay System	100 assays	E1910
Dual-Luciferase® Reporter Assay System 10-Pack	10 x 100 assays	E1960
Dual-Luciferase® Reporter Assay System	1000 assays	E1980
Dual-Glo® Luciferase Assay System	10 ml	E2920
Dual-Glo® Luciferase Assay System	100 ml	E2940
Dual-Glo® Luciferase Assay System	10 x 100 ml	E2980

转染试剂

产品	规格	目录号
TransFast™ Transfection Reagent	1.2 mg	E2431
FuGENE® HD Transfection Reagent	1 ml	E2311
FuGENE® HD Transfection Reagent	5 x 1 ml	E2312
FuGENE® 6 Transfection Reagent	1 ml	E2691
FuGENE® 6 Transfection Reagent	5 x 1 ml	E2692
FuGENE® 6 Transfection Reagent	0.5 ml	E2693
ViaFect™ Transfection Reagent	0.75 ml	E4981
ViaFect™ Transfection Reagent	2 x 0.75 ml	E4982

www.promega.com.cn/products/luciferase-assays/



关注 Promega 微信公众号



价格查询



中文说明书



实验工具



技术资料



市场活动



经销商信息

普洛麦格 (北京) 生物技术有限公司

Promega (Beijing) Biotech Co., Ltd

地址: 北京市东城区北三环东路 36 号环球贸易中心 B 座 907-909

电话: 010-58256268

网址: www.promega.com

技术支持电话: 400 810 8133

技术支持邮箱: chinatechserv@promega.com

更新时间: 2022.3