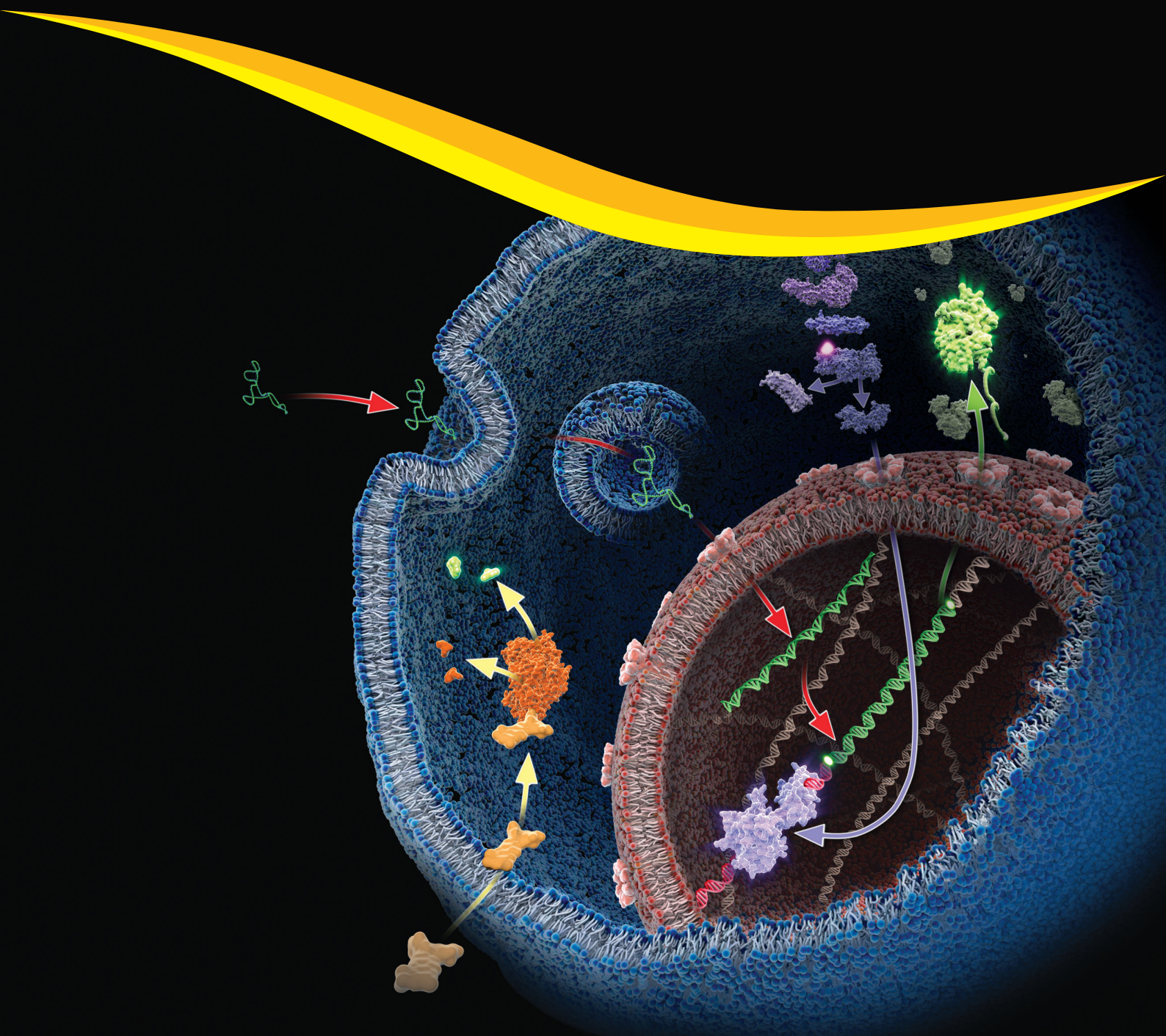


miRNA 及 siRNA 功能研究技术

pmirGLO 萤光素酶报告基因技术

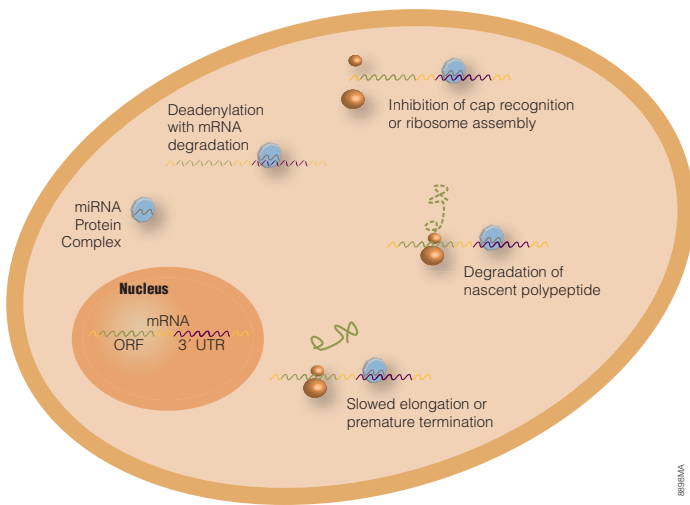


微小 RNA

microRNA(miRNA)

microRNA(miRNA) 即微小 RNA 广泛存在于真核细胞中, 是内源性的具有调控功能的非编码 RNA(长度约 17~24 个核苷酸), 已知其在动植物中介导广泛的生物学功能。在人类中, miRNA 参与发育、分化及代谢等多方面的调控, 并被证实调控从神经元发育到干细胞分裂等多种过程的基因表达, 并且 miRNA 表达失调与多种疾病状态相关, 包括神经退行性疾病、心血管疾病和癌症。但到目前为止, 只有一小部分 miRNA 生物学功能得到阐明。更多的 miRNA 功能还尚待研究。

通常一个 miRNA 分子可以靶向多个靶 mRNA, 一个靶 mRNA 可以被多个不同 miRNA 协同调节, 只结合一个 miRNA 无法充分发挥抑制翻译或诱导 mRNA 降解的调节作用, 通常需要多个 miRNA 结合至同一靶 mRNA。由于 miRNA 作用的多样性和复杂性, 造成了对 miRNA 功能研究平台有更高的要求。



miRNA 的作用原理及报告基因的应用

miRNA 能够形成 RNA 诱导的沉默复合体 (RNA-induced silencing complex, RISC), 主要通过碱基互补配对的方式结合靶 mRNA 的 3'非翻译区 (UTR) 的与种子序列互补的片段 (miRNA 中介导靶序列结合的 2-7 个核苷酸组成的小片段区域称为 “seed sequence”, 即种子序列), 根据互补程度的不同降解靶 mRNA 或者阻遏靶 mRNA 的翻译, 使靶基因沉默。

研究 miRNA 调控作用的实验常采用基于细胞的报告基因系统, 萤光素酶 (萤光素酶) 就是最为经典的报告基因技术之一。通常实验中, 报告基因被克隆至待研究的 3'-UTR 序列上游, 并与 miRNA 共转染至培养细胞中。

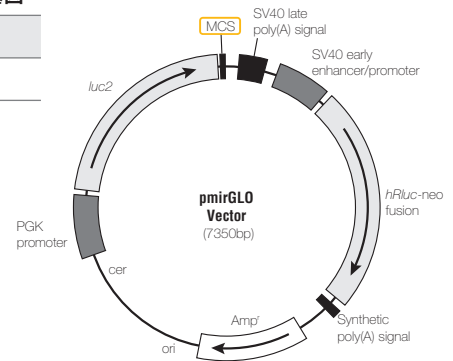
miRNA 功能研究思路举例

- 1 预测结合靶基因的可能 miRNA (通过预测网站进行预测)
- 2 RNA 或蛋白检测验证靶基因与 miRNA 表达水平存在负相关
- 3 通过报告基因在细胞水平确认位点

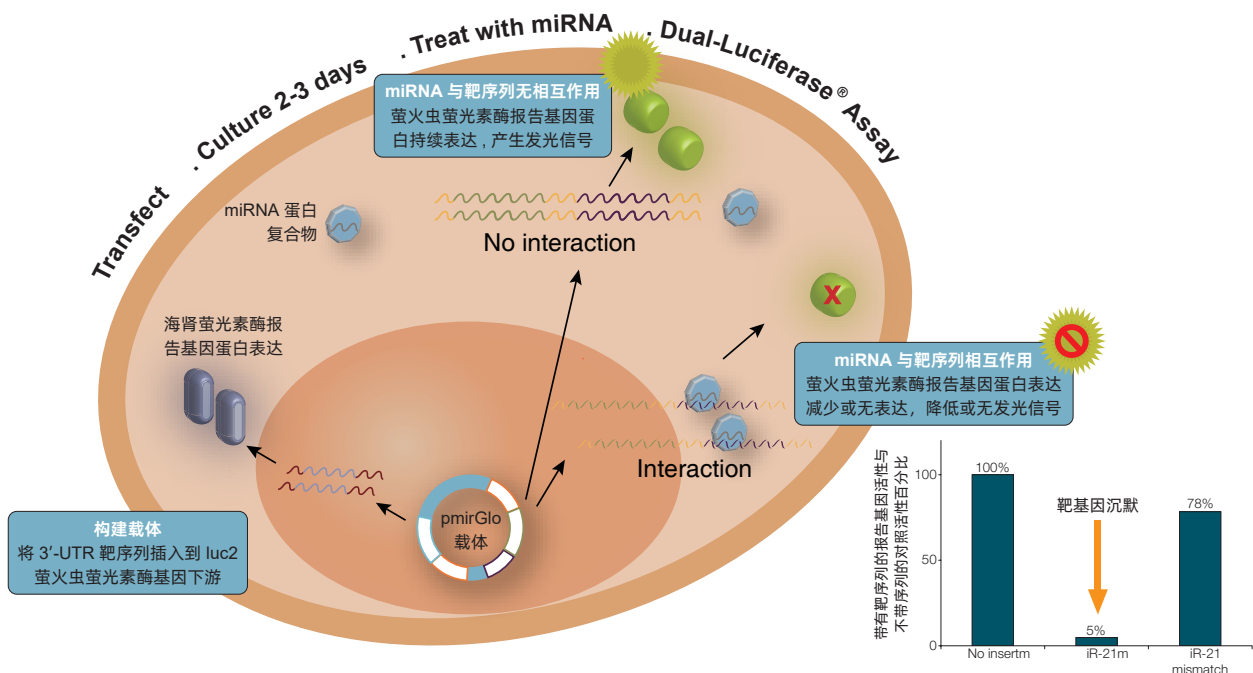
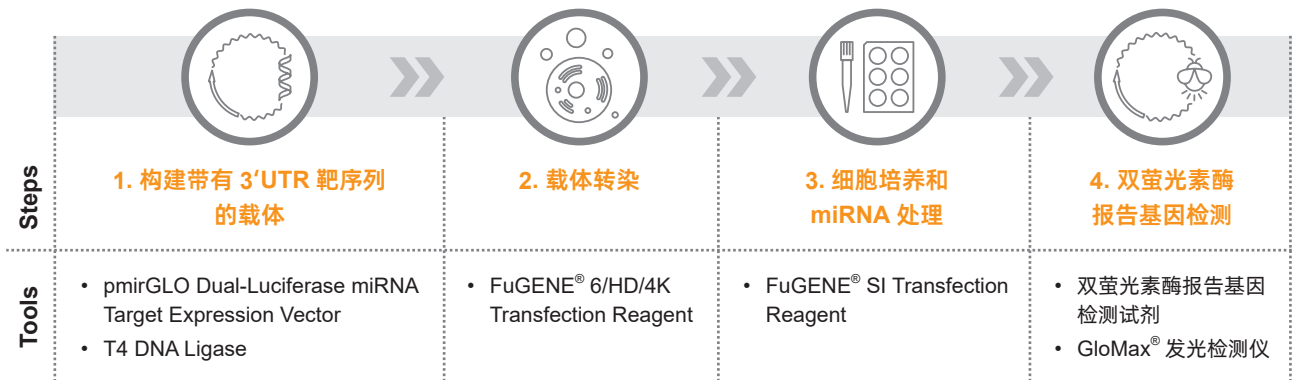
pmirGLO 报告基因技术

pmirGLO 报告基因检测技术是由 pmirGLO 载体，转染试剂和检测试剂及配套相关产品组成的完整解决方案。pmirGlo 报告基因载体是在一个载体上带有两个报告基因；一个是由组成型 PGK 启动子驱动的主报告基因，研究者可将 miRNA 结合的 3'-UTR 靶序列插入到主报告基因下游或 3'端，而另外一个内参报告基因，由组成型 SV40 启动子驱动表达。如 miRNA 与靶序列结合发挥作用，则主报告基因表达会下降，而内参报告基因不受影响。Promega 提供 2 种报告基因组合的载体产品，供您挑选。

载体	主报告基因	内参报告基因
pmirGLO Dual-Luciferase miRNA Target Expression Vector	Firefly	Renilla
pmirGLO NanoGlo Vector	NanoLuc	Firefly



- **细胞水平的验证方法：**更能说明 miRNA 对其靶序列或靶基因的作用。
- **久经验证的报告基因检测法：**拥有大量应用文献。
- **适用于内源 miRNA 活性研究：**主报告基因上游带有中等转录活性的 PGK 启动子，对基因表达抑制反应更灵敏，可用于内源性 miRNA 活性的检测。
- **携带新霉素抗性：**可用于建立稳转细胞株。



上图：使用 pmirGLO Dual-Luciferase miRNA 载体进行实验的操作流程示例图

产品列表

产品应用	产品名称	规格	目录号
用于 3'UTR 报告 基因载体克隆	pmirGLO Dual-Luciferase miRNA Target Expression Vector	20 ug	E1330
	pmirGLO NanoGlo miRNA Target Expression Vector	20 ug	请咨询 Promega
	T4 DNA Ligase	100 u	M1801
		500 u	M1804
T4 DNA Ligase(HC)	500 u	M1794	
	PureYield™ Plasmid Miniprep System	250 prep	A1222
		100 prep	A1223
载体转染	FuGENE® 6 Transfection Reagent	1 ml	E2691
		5 × 1 ml	E2692
		0.5 ml	E2693
	FuGENE® HD Transfection Reagent	1 ml	E2311
		5 × 1 ml	E2312
	FuGENE® 4K Transfection Reagent	1 ml	E5911
5 × 1 ml		E5912	
miRNA 转染	FuGENE® SI Transfection Reagent	1 ml	E9311
		5 × 1 ml	E9312
		0.2 ml	E9321
转染毒性检测 *	CellTiter-Glo® Luminescent Cell Viability Assay	10 ml	G7570
		10 × 10 ml	G7571
		100 ml	G7572
		10 × 100 ml	G7573
双萤光素酶 报告检测试剂 (Firefly/Renilla)	Dual-Luciferase® Reporter Assay System	100 assays	E1910
		10 × 100 assays	E1960
		1000 assays	E1980
	Dual-Glo® Luciferase Assay System	10 ml	E2920
		100 ml	E2940
双萤光素酶 报告检测试剂 (NanoLuc/Firefly)	Nano-Glo® Dual-Luciferase® Reporter Assay System	10 ml	N1610
		100 ml	N1620
		10 × 10 ml	N1630
检测仪器	GloMax® Discover System	1 台	GM3000
其他相关产品			
反转录	GoScript™ Reverse Transcription System	50 rxns	A5000
		100 rxns	A5001
	Eastep™ RT Master Mix Kit	50 rxns	LS2050
		200 rxns	LS2052
Eastep™ RT Master Mix Kit, with No-RT Control	200 rxns	LS2054	
qPCR	GoTaq® qPCR Master Mix	1 ml	A6000
		5 ml (500 × 20 ul rxns)	A6001
		25 ml (2500 × 20 ul rxns)	A6002
	Eastep™ qPCR Master Mix Kit	500 rxns	LS2062
2000 rxns		LS2068	

* 除了 CellTiter-Glo 试剂，Promega 还可提供多种细胞活力和增殖检测试剂，详情请咨询 Promega

应用文献汇总

应用方向	文章信息	应用产品
miRNA 与疾病机制研究	<p>OncomiR-181a promotes carcinogenesis by repressing the extracellular matrix proteoglycan decorin in hepatocellular carcinoma. BMC Gastroenterol. 2024 Sep 30;24(1):337. DOI: 10.1186/s12876-024-03413-6.</p>	<ul style="list-style-type: none"> • pmirGLO Dual-Luciferase miRNA Target Expression Vector • Steady-Glo® Luciferase Assay System
	<p>Hsa-miR-874-3p Reduces Endogenous Expression of RGS4-1 Isoform In Vitro. Genes (Basel). 2024 Aug 11;15(8):1057. DOI: 10.3390/genes15081057.</p>	<ul style="list-style-type: none"> • pmirGLO Dual-Luciferase miRNA Target Expression Vector • Dual-Luciferase® Reporter Assay System
	<p>miR-22 and miR-29a Are Members of the Androgen Receptor Cistrome Modulating LAMC1 and Mcl-1 in Prostate Cancer Mol Endocrinol. 2015 Jun 8;29(7):1037–1054. DOI: 10.1210/me.2014-1358</p>	<ul style="list-style-type: none"> • pmirGLO NanoGlo miRNA Target Expression Vector • Nano-Glo® Dual-Luciferase® Reporter Assay System • anti-PARP p85
miRNA 调控网络与相互作用研究	<p>miR-122-5p is involved in posttranscriptional regulation of the mitochondrial thiamin pyrophosphate transporter (<i>SLC25A19</i>) in pancreatic acinar cells. Am J Physiol Gastrointest Liver Physiol. 2023 Aug 2;325(4):G347–G355. DOI: 10.1152/ajpgi.00106.2023</p>	<ul style="list-style-type: none"> • pGEM-T easy vector • pmirGLO Dual-Luciferase miRNA Target Expression Vector • Dual-Luciferase® Reporter Assay System
	<p>Mouse Pxt1 expression is regulated by Mir6996 miRNA. Theriogenology. 2023 Oct 15;210:9-16. DOI: 10.1016/j.theriogenology.2023.07.010</p>	<ul style="list-style-type: none"> • pGEM-T Easy Vector • pmirGLO Dual-Luciferase miRNA Target Expression Vector • Dual-Glo® Luciferase Assay System
	<p>miR-29a, b, and c regulate SLC5A8 expression in intestinal epithelial cells. Am J Physiol Gastrointest Liver Physiol. 2021 Aug 1;321(2):G223-G231. DOI: 10.1152/ajpgi.00148.2021</p>	<ul style="list-style-type: none"> • pmirGLO Dual-Luciferase miRNA Target Expression Vector • Dual-Glo® Luciferase Assay System • GloMax 2020 Luminometer
miRNA 靶向治疗开发	<p>MicroRNA-205-5p: A potential therapeutic target for influenza A. J Cell Mol Med. 2022 Dec;26(23):5917-5928. DOI: 10.1111/jcmm.17615</p>	<ul style="list-style-type: none"> • pmirGLO Dual-Luciferase miRNA Target Expression Vector • Dual-Luciferase® Reporter Assay System
	<p>Downregulation of NHE-3 (SLC9A3) expression by MicroRNAs in intestinal epithelial cells. Am J Physiol Cell Physiol. 2022 Oct 3;323(6):C1720–C1727. doi: 10.1152/ajpcell.00294.2022</p>	<ul style="list-style-type: none"> • pmirGLO Dual-Luciferase miRNA Target Expression Vector • Dual-Luciferase® Reporter Assay System • (HRP)-conjugated goat anti-rabbit IgG antibody
miRNA 作为诊断与预后标志物研究	<p>DLEU7-AS1 promotes renal cell cancer by silencing the miR-26a-5p/coronin-3 axis. Clin Kidney J. 2022 Feb 28;15(8):1542-1552. DOI: 10.1093/ckj/sfac061</p>	<ul style="list-style-type: none"> • pmirGLO Dual-Luciferase miRNA Target Expression Vector • Dual-Glo® Luciferase Assay System
疫苗研发	<p>Implementation of immuno-chemoinformatics approaches to construct multi-epitope for vaccine development against Omicron and Delta SARS-CoV-2 variants Vacun. 2022 Sep-Dec;23:S18-S31. DOI: 10.1016/j.vacun.2022.05.006</p>	<ul style="list-style-type: none"> • pmirGLO Dual-Luciferase miRNA Target Expression Vector
疾病机制研究	<p>Roles of phosphatidylinositol-3-kinases signaling pathway in inflammation-related cancer: Impact of rs10889677 variant and buparlisib in colitis-associated cancer. World J Gastroenterol. 2023 Oct 28;29(40):5543-5556. DOI: 10.3748/wjg.v29.i40.5543.</p>	<ul style="list-style-type: none"> • GoTaq® Green Master Mix • PureYield™ Plasmid Miniprep System • pmirGLO Dual-Luciferase miRNA Target Expression Vector • Dual-Glo® Luciferase Assay System • GloMax 2020 Luminometer

miRNA 及 siRNA 功能研究技术



关注 Promega 生命科学公众号，您可获得



产品信息



价格查询



中文说明书



讲座视频



技术资料



实验工具



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：2025.04