



HiBiT 蛋白标签系统使用指南:

Step1. 如何用 HiBiT 标签标记蛋白

方法 1: 构建载体进行重组表达			方法 2: 用 CRISPR/Cas9 导入 HiBiT 内源标签		
• 可选载体: - HiBiT CMV-neo Flexi[®] Vectors 载体系列 - pBiT3.1 HiBiT MCS Vectors 载体系列			① 选择即用型细胞系 Promega 提供 100+ 针对热门靶标的即用型细胞系。  扫码查看完整细胞系列表	② 定制细胞系 联系 Promega 定制满足您研究需求的细胞系。 https://www.promega.com.cn/custom-solutions/tailored-solutions/	③ 自行 DIY 参考 Promega 提供的操作指导自行完成 CRISPR 基因敲入。 https://www.promega.com.cn/applications/crispr-diy-overview/
• 可选转染试剂					
FugENE[®] HD E2311 1ml E2312 5 x 1ml	FugENE[®] 6 E2691 1ml E2692 5 x 1ml E2693 0.5ml	ViaFect[™] E4981 0.75ml E4982 2 x 0.75ml E4983 0.2ml			

Step 2. 如何检测 HiBiT 标签标记的蛋白

 裂解型检测	 细胞外检测	 细胞内检测	 印迹检测	 基于抗体的检测
裂解检测试剂定量 HiBiT 标签蛋白的总量。	定量活细胞表面或分泌型表达的蛋白。	在细胞内共表达 LgBiT, 定量细胞内的 HiBiT 标签蛋白; 无需裂解细胞; 实时动力学研究可达 72 小时。	使用简单的发光信号即可在几分钟内检测印迹膜上任意标签化蛋白。	执行传统的基于抗体的检测, 例如通过 WB, IP, ICC, pull-down 的方法检测 HiBiT 标签蛋白。
N3030 10ml N3040 100ml N3050 10 x 100ml	N2420 10ml N2421 100ml N2422 10 x 100ml	Nano-Glo[®] 活细胞检测系统 N2011 100 assays 长时程动态监测底物 N2570 0.1ml N2580 0.1ml N2590 0.2ml LgBiT mRNA 递送系统 NE1120 10plates LgBiT 表达载体 N2681 20µg LgBiT 稳转细胞系 N2672 1each	N2410 100ml	Anti-HiBiT 单抗 N7200 100µg N7210 5 x 100µg Anti-HiBiT 单抗, FarRed647 N7221 50µg Anti-HiBiT 单抗, Green488 N7231 50µg Anti-HiBiT Magne[®] 磁珠 N7300 1mL N7301 5mL N7400 0.1mL

注: 以上列出的产品均可通过目录号前往 www.promega.com.cn 或 ministe.promega.com.cn 查询产品信息及更多相关资源。



HiBiT 蛋白标签高分引文列表

应用方向	引用文献	主要使用产品
靶蛋白降解	Targeted protein degradation via intramolecular bivalent glues. Hsia, O., <i>et al.</i> (2024) <i>Nature</i> 627, 204–211. https://doi.org/10.1038/s41586-024-07089-6	Nano-Glo® HiBiT Lytic detection system (Cat# N3030) GloMax® Discover microplate reader (Cat# GM3000)
生物发光成像	An improved pathway for autonomous bioluminescence imaging in eukaryotes. Shakhova, E.S., <i>et al.</i> (2024) <i>Nat Methods</i> 21, 406–410. https://doi.org/10.1038/s41592-023-02152-y	Nano-Glo® HiBiT Lytic detection system (Cat# N3030)
GPCR	An inverse agonist of orphan receptor GPR61 acts by a G protein-competitive allosteric mechanism. Lees, J.A., <i>et al.</i> (2023) <i>Nat Commun</i> 14, 5938. https://doi.org/10.1038/s41467-023-41646-3	Nano-Glo® HiBiT Extracellular Detection System (Cat# 2421) Nano-Glo® HiBiT Lytic detection system (Cat# N3030)
GPCR	Structural insights into constitutive activity of 5-HT6 receptor. Licong He, <i>et al.</i> (2023) <i>PNAS</i> 120 (14) , e2209917120. https://doi.org/10.1073/pnas.2209917120	Nano-Glo® HiBiT Extracellular Detection System (Cat# 2421)
认知障碍炎症	Platelet factors attenuate inflammation and rescue cognition in ageing. Schroer, A.B., <i>et al.</i> (2023) <i>Nature</i> 620, 1071–1079. https://doi.org/10.1038/s41586-023-06436-3	Nano-Glo® HiBiT Lytic detection system (Cat# N3030)
蛋白递送基因治疗	Programmable protein delivery with a bacterial contractile injection system. Kreitz, J., <i>et al.</i> (2023) <i>Nature</i> 616, 357–364. https://doi.org/10.1038/s41586-023-05870-7	Nano-Glo® HiBiT blotting system (Cat# N2410)
药物研发炎症检测	Probing expression of E-selectin using CRISPR-Cas9-mediated tagging with HiBiT in human endothelial cells. Lydia Ogrodzinski, <i>et al.</i> (2023) <i>iScience</i> 26(7):107232. doi: 10.1016/j.isci.2023.107232 .	Nano-Glo® HiBiT Lytic Detection System (Cat# N3040) Anti-HiBiT Monoclonal Antibody (Cat# N7200)
病毒研究	Virological characteristics of the SARS-CoV-2 Omicron BA.2 spike Yamasoba, Daichi <i>et al.</i> (2022) <i>Cell</i> 185(12), 2103 - 2115.e19 DOI:https://doi.org/10.1016/j.cell.2022.04.035	Nano-Glo® HiBiT lytic detection system (Cat# N3040) GloMax® Explorer Microplate Reader (Cat# GM3510)
靶蛋白降解	Antibody targeting of E3 ubiquitin ligases for receptor degradation. Marei, H., <i>et al.</i> (2022) <i>Nature</i> 610, 182–189. https://doi.org/10.1038/s41586-022-05235-6	Nano-Glo® HiBiT Extracellular Detection System (Cat# 2421)
糖酵解荧光成像	Hexokinase 3 enhances myeloid cell survival via non-glycolytic functions. Seiler, K., <i>et al.</i> (2022) <i>Cell Death Dis</i> 13, 448. https://doi.org/10.1038/s41419-022-04891-w	Anti-HiBiT Monoclonal Antibody (Cat# N7200)
药物研发	Determining ADCC activity of antibody-based therapeutic molecules using two bioluminescent reporter-based Bioassays. Garvin, D. <i>et al.</i> (2021) <i>Curr Protoc.</i> 1(11):e296 https://doi.org/10.1002/cpz1.296	Nano-Glo® HiBiT Extracellular Detection System (cat# N2420; N2421; N2422)

普洛麦格 (北京) 生物技术有限公司
Promega (Beijing) Biotech Co., Ltd

技术服务电话: 400 810 8133

技术服务邮箱: chinatechserv@promega.com

官网: www.promega.com 微网站: wechat.promega.com.cn

邮箱: marketing@promega.com.cn

更新时间: 2026.07

关注 Promega
生命科学



联系
授权经销商

