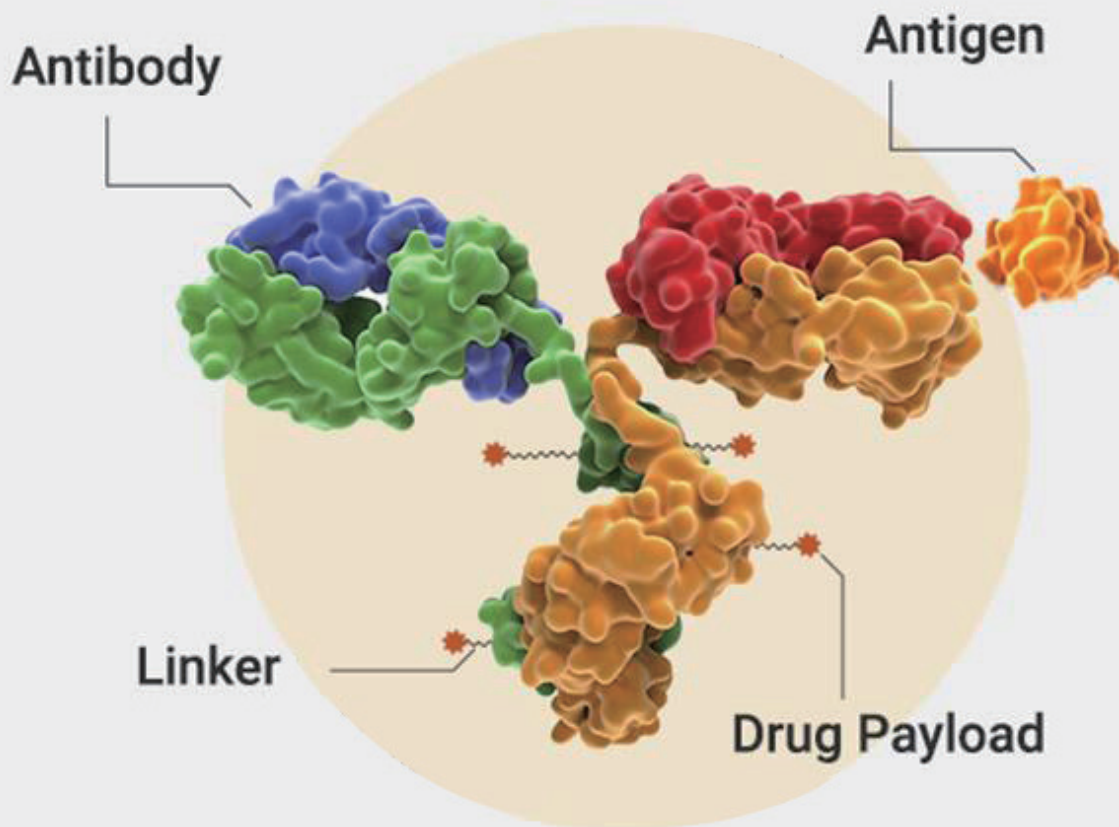


ADC 药物研发解决方案

抗体筛选及表征 | Payload 功能分析 | 抗体偶联技术开发
Fc 受体表征 | 抗体内化 | 靶细胞杀伤 | 细胞因子检测



ADC 药物研发解决方案矩阵

1. 药物发现

支持您在选择合适的靶点、抗体、连接器 (Linker) 和小分子细胞毒药物 (Payload), 以及开发有效的偶联技术等方面进行高效研究。

2. 工艺开发

多种抗体表征分析工具组合, 帮助您在 ADC 药物研发中进行严格的质量控制检测。

3. 临床前研究

Promega 拥有丰富的抗体药物研发用工具, 帮助您在临床前研究中全方面评估 ADC 药物的安全性和有效性。

Payload 功能分析



细胞毒性和凋亡检测系统

精确评估有效载荷诱导的细胞毒性效应和凋亡, 指导开发更有效的癌症治疗方法。



细胞健康
解决方案



免疫原性细胞死亡 (ICD) 检测系统

Lumit[®] 免疫检测法能够对标志着 ICD 的关键细胞因子进行快速、精确可靠的定量。



Lumit[®] 技术
解决方案



降解剂分析检测技术

基于活细胞研究的解决方案, 用于分析降解剂或 PROTACs 作为 ADC 有效载荷的功能, 以便仔细评估降解剂有效载荷的效力和作用机制。



蛋白降解
解决方案



寡核苷酸分析工具

靶向 RNA 研究工具能够详细分析针对 RNA 的有效载荷作用及其对药物开发的影响。



靶向 RNA
研究工具

ADC 开发及表征

抗体表征工具可帮助确认生物治疗性蛋白质序列, 包括 ADC 的偶联位点, 评估 ADC 的稳定性和降解情况。揭示药物与抗体的比例。

Magne[®] Protein G/A 磁珠使偶联和纯化过程变得直接、可扩展且快速。允许在珠上直接进行偶联。



抗体表征工具

质量控制



质谱分析

- 监测宿主细胞蛋白杂质;
- 确保每批 ADC 都符合严格的质量标准及批次间的一致性。



抗体表征
工具



抗体内化分析

Antibody Internalization Bioassay 的数据结果可准确反映抗原表达与 ADC 结合动力学, 定量并监测抗体内化, 以筛选最优抗体设计。



抗体内化
新品介绍

pH 敏感性染料能够精确追踪抗体进入细胞的内吞过程, 有助于评估和增强 ADC 的治疗潜力。



pHAb dye
产品介绍

药效与机制研究



Fc 受体表征工具

Fc 受体 (FcR) 的相互作用会直接影响 ADC 药物作用机制及免疫系统的参与。针对 ADCC、ADCP 以及 CDC 的检测工具, 可用于优化 ADC 的疗效。



Fc 受体功能
分析与表征



细胞杀伤检测系统

HiBiT 靶细胞杀伤生物测定法能够在混合培养中观察特定类型细胞的死亡。有助于分析 ADC 分子的旁观者杀伤效应和特异性。



HiBiT 蛋白
标签技术



活体成像检测技术

通过活体成像实验, 监测肿瘤的生长和转移, 以此评估 ADC 药物的疗效和了解肿瘤的生物行为。



小动物活体
成像解决方案

ADC 药物研发工具列表

应用方向	主要优势 / 特点	产品名称	规格	目录号
抗体内化检测	- 免洗涤、快速定量 - 抗原特异性内化 - 检测不受 Payload 干扰 - 广泛的适用性 (贴壁细胞、悬浮细胞)	Antibody Internalization Bioassay 	1 each	CS372004
		Antibody Internalization Bioassay 10X	1 each	CS372008
	- pH 感应式染料标记抗体 - 加样读数、无洗涤步骤、操作简单 - 信噪比更高 - 实时动态监测	pHAb Amine Reactive Dye	1 × 250 µg	G9841
		pHAb Thiol Reactive Dye	1 × 250 µg	G9831
抗体偶联技术开发	用于抗体捕获的高特异性、高容量亲和磁珠	Magne™ Protein G Beads	1 ml	G7471
		Magne™ Protein A Beads	1 ml	G8781
细胞活力检测	- 金标准检测法 - 快速, 最快 10 分钟孵育 - 灵敏, 可检测低至 10 个细胞 - 高通量, 轻松兼容 96-1536 通量	CellTiter-Glo® Luminescent Cell Viability Assay(CTG)	10 ml	G7570
	- CTG 的迭代产品 - 单一试剂, 可直接使用 - 试剂稳定时间更长	CellTiter-Glo® 2.0 Assay	10 ml	G9241
	- 专为 3D 培养样本而设计 - 具有 CTG 产品所有特点	CellTiter-Glo® 3D Cell Viability Assay	10 ml	G9681
细胞凋亡检测	- 适用于多种类型培养细胞 - 实时检测剂量依赖和时间依赖的凋亡进程 - 可与其他检测试剂叠加使用	RealTime-Glo™ Annexin V Apoptosis and Necrosis Assay	100 assays	JA1011
免疫原性细胞死亡: Lumit® 细胞因子检测	- 简单的“加样 - 读数”模式, 无需洗涤, 无需固定 - 操作最少只需 70min - 生物发光法, 仅需发光检测仪 - 适用于 96 或 384 孔板高通量检测	Lumit® IL-2 (Human) Immunoassay	100 assays	W6020
			1,000 assays	W6021
			5 x 100 assays	W6022
		Lumit® IL-6 (Human) Immunoassay	100 assays	W6030
			1,000 assays	W6031
			5 x 100 assays	W6032
		Lumit® IL-10 (Human) Immunoassay	100 assays	W6070
			1,000 assays	W6071
			5 x 100 assays	W6072
		Lumit® IFN-γ (Human) Immunoassay	100 assays	W6040
			1,000 assays	W6041
			5 x 100 assays	W6042

应用方向	主要优势 / 特点	产品名称	规格	目录号
免疫原性细胞死亡： Lumit® 免疫检测	- 简单的“加样 - 读数”模式，无需洗涤，无需固定 - 操作最少只需 70min - 生物发光法，仅需发光检测仪 - 适用于 96 或 384 孔板高通量检测	Lumit® TNF-α (Human) Immunoassay	100 assays	W6050
			1,000 assays	W6051
			5 x 100 assays	W6052
		Lumit® IL-4 (Human) Immunoassay	100 assays	W6060
			1,000 assays	W6061
			5 x 100 assays	W6062
		Lumit® Human IL-1β Immunoassay	100 assays	W6010
			1,000 assays	W6011
			5 x 100 assays	W6012
		Lumit® Mouse IL-1β Immunoassay	100 assays	W7010
			1,000 assays	W7011
			5 x 100 assays	W7012
		Lumit® IL-8 (Human) Immunoassay	100 assays	W1460
			500 assays	W1462
			1,000 assays	W1461
		Lumit® Active IL-18 (Human) Immunoassay	100 assays	W1910
			1,000 assays	W1911
			5 x 100 assays	W1912
		Lumit® Human IL-18 Standard	25 µl	W191A-C
		Lumit® IFN-β (Human) Immunoassay	100 assays	W1810
			1,000 assays	W1811
			5x 100 assays	W1812
		Human IFN-β Standard	25 µl	W149A-C
		Lumit® IL-17A (Human) Immunoassay	100 assays	W1430
			1,000 assays	W1431
			5 x 100 assays	W1432
		Human IL-17A Standard	25 µl	W143A-C
		Lumit® IL-12 p70 (Human) Immunoassay	100 assays	W1850
			500 assays	W1852
			1,000 assays	W1851
			25 µl	W185A-C

应用方向	主要优势 / 特点	产品名称	规格	目录号
旁观者效应： 靶细胞杀伤检测	- 可以终点法或动态检测 - 能够进行长达 24 小时或更长时间的分析 - 灵敏、简单，可用于研发和 / 或 QC Lot Release	H929 (HiBiT) TCK Bioassay, Propagation Model	2 vials	M2412
		Raji (HT-HiBiT) TCK Bioassay, Propagation Model	2 vials	M2452
		Ramos (HiBiT) TCK Bioassay, Propagation Model	2 vials	M2472
		其他 HiBiT TCK Bioassays (如 A549 cells, SK-BR-3 cells, U937 cells, A375 cells, T2 cells 等等)	Please enquire	
Fc 受体表征： ADCC 活性检测	- 基于 MoA 的报告基因检测法 - ICH 指南预验证 - 操作简单，可高通量操作 - 误差小，准确性好 - 可用于批签发	ADCC Bioassay Effector Cells, Propagation Model	1 each	G7102
Fc 受体表征： ADCP 活性检测	- 基于 MoA 的报告基因检测法 - ICH 指南预验证 - 操作简单，可高通量操作 - 误差小，准确性好 - 可用于批签发	ADCP Reporter Bioassay (THP-1)	1 each	GA1272
Fc 受体表征： CDC 活性检测	- 可反映细胞膜完整性 10 个死细胞即可检测 15 分钟孵育即可	CytoTox-Glo™ Cytotoxicity Assay	10ml	G9290
	- 评估 C1q 抗体结合以评估补体依赖性 - ELISA 替代方案，免洗，快速，无需固定，兼容高通量 60 分钟内即可获得结果	Lumit® C1q Binding Assay 	1 each	CS3695A04
Fc 受体表征： Fc 受体结合检测	- 所需样本量少 (10-25 µl) - 操作简单，无需洗涤和转移 1 小时内完成检测 - 兼容高通量	FcRn Binding Immunoassay	100 assays	W1151
		FcRn Binding Immunoassay 10X	1,000 assays	W1152
		FcγRI Binding Immunoassay	100 assays	W7080
		FcγRI Binding Immunoassay 10X	1,000 assays	W7081
		FcγRIIa (H131) Binding Immunoassay	100 assays	W7070
		FcγRIIa (H131) Binding Immunoassay 10X	1,000 assays	W7071
		FcγRIIa (R131) Binding Immunoassay	100 assays	W7060
		FcγRIIa (R131) Binding Immunoassay 10X	1,000 assays	W7061

应用方向	主要优势 / 特点	产品名称	规格	目录号
Fc 受体表征： Fc 受体结合检测	- 所需样本量少（10-25 µl） - 操作简单，无需洗涤和转移 1 小时内完成检测 - 兼容高通量	FcγRIIIa (V158) Binding Immunoassay	100 assays	W7050
		FcγRIIIa (V158) Binding Immunoassay 10X	1,000 assays	W7051
		FcγRIIIa (F158) Binding Immunoassay	100 assays	W7040
		FcγRIIIa (F158) Binding Immunoassay 10X	1,000 assays	W7041
		FcγRIIb Binding Immunoassay	100 assays	W7030
		FcγRIIb Binding Immunoassay 10X	1,000 assays	W7031
质谱分析	用于肽图谱和蛋白质组学分析的重组胰凝乳蛋白酶	rChymoSelect™ MS Grade Protease Kit 	25 µg	CS333204
			100 µg	CS333205
	高度特异性免疫球蛋白 IgG 降解酶	IdeS/IdeZ Protease	5,000 units	V7511/ V8341
	重组糖苷酶	PNGase F	500 u	V4831
	胰蛋白酶	Sequencing Grade Modified Trypsin	5 × 20 µg	V5111
	质谱级重组胰蛋白酶	Trypsin Platinum, Mass Spectrometry Grade	100 µg	VA9000
	Trypsin 与 Lys-C 的混合物	Trypsin/Lys-C Mix, Mass Spec Grade	20 µg	V5071
	特异性替代蛋白酶	ProAlaase, Mass Spec Grade	5 µg	VA2161
		Asp-N, Sequencing Grade	2 µg	V1621
		Glu-C, Sequencing Grade	5 × 10 µg	V1651
	低特异性替代蛋白酶	Chymotrypsin, Sequencing Grade	25 µg	V1061
	非特异性蛋白酶	Pepsin	250 mg	V1959
	低 PH 蛋白酶	AccuMAP™ Low pH Protein Digestion Mini Kit	10 reactions	VA1040

注：列表中的产品均有其他规格提供，请浏览 [wechat.Promega.com.cn](https://wechat.promega.com.cn) 查看更多详情。

ADC 药物研发应用文献列表

应用方向	文章信息	应用产品
BET 蛋白降解剂， 肿瘤研究	Nakazawa, Y. <i>et al.</i> (2024) Delivery of a BET protein degrader via a CEACAM6-targeted antibody–drug conjugate inhibits tumour growth in pancreatic cancer models. Nature.Communications . 15:2192. DOI:10.1038/s41467-024-46167-1 .	• Nano-Glo® Extended Live Cell Substrates • CellTiter-Glo® 3D Cell Viability Assay • VivoGlo™ Luciferin, In Vivo Grade • Dual-Glo® Luciferase Assay System • Maxwell® RSC simplyRNA Tissue Kit
ADC 筛选工具开发	Vasic, V., <i>et al.</i> (2024) Generation of binder-format-payload conjugate-matrices by antibody chain-exchange. Nat Commun 15, 9406. https://doi.org/10.1038/s41467-024-53730-3	• pHAb Amine Reactive Dye Kit
DNA 损伤修复， 白血病研究	Ghelli Luserna di Rorà, A., <i>et al.</i> (2024) Exploring the role of PARP1 inhibition in enhancing antibody–drug conjugate therapy for acute leukemias: insights from DNA damage response pathway interactions. J Transl Med 22, 1062. https://doi.org/10.1186/s12967-024-05838-9	• RealTime-Glo® MT Cell Viability Assay
霍奇金淋巴瘤 (HL)	F Morelli, <i>et al.</i> (2024) Antibody-Drug Conjugate Made of Zoledronic Acid and the Anti-CD30 Brentuximab-Vedotin Exert Anti-Lymphoma and Immunostimulating Effects. Cells 13(10):862. https://doi.org/10.3390/cells13100862	• CellTiter-Glo® Luminescent Cell Viability Kit • Caspase Glo®-3/7 3D Assays • Caspase-Glo® 9 Assay Systems
GPCR 与结直肠癌研究	Jacob, J., <i>et al.</i> (2023). An antibody-drug conjugate targeting GPR56 demonstrates efficacy in preclinical models of colorectal cancer. British journal of cancer , 128(8), 1592–1602. https://doi.org/10.1038/s41416-023-02192-3	• Reporter vector pGL4.34[luc2P/SRF-RE/Hygro] • CellTiter-Glo® assay
癌症免疫治疗	Sue M, <i>et al.</i> (2024) Blockade of SIRPα-CD47 axis by anti-SIRPα antibody enhances anti-tumor activity of DXd antibody-drug conjugates. PLoS One . Jun 6;19(6):e0304985. https://doi.org/10.1371/journal.pone.0304985	• FuGENE® 6 transfection reagent • Lumit® HMGB1 Immunoassay • RealTime-Glo Extracellular ATP Assay • BacTiter-Glo Reagent • CellTiter-Glo® Luminescent Cell Viability Assay

注：以上应用产品可前往 www.promega.com.cn 或者 wechat.promega.com.cn 查询更多信息。

ADC 药物研发

www.promega.com/applications/biologics-drug-discovery/antibody-drug-conjugate/



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：2026.09