

Double-stranded RNA (dsRNA) ELISA Kit (J2 based)

Cat.No. L01020

产品特点



双J2抗体夹心ELISA法

- J2 抗体既是捕获抗体，又是检测抗体



快速检测时间

- 3 hours



宽线性范围

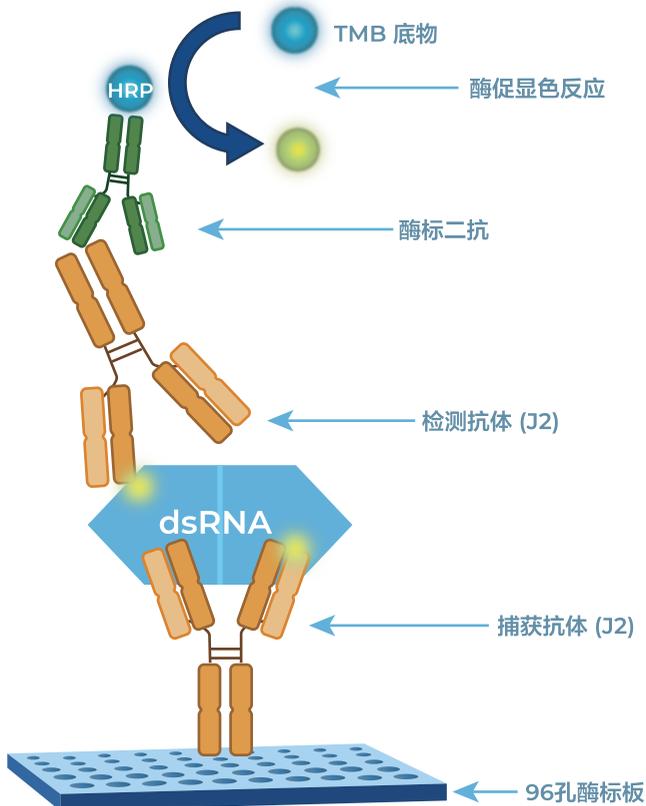
- 0.123-30 ng/mL



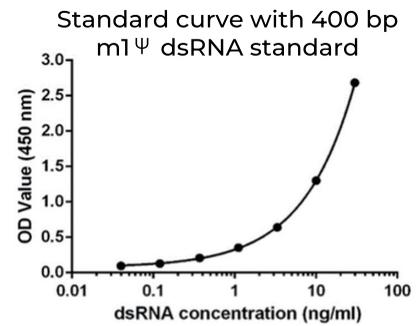
标准品

- 400 bp 1mΨ修饰的dsRNA
- 金斯瑞可以提供不同长度和不同修饰的标准品
- *标准品参考Moderna发表文献中的长度和工艺进行合成

检测原理

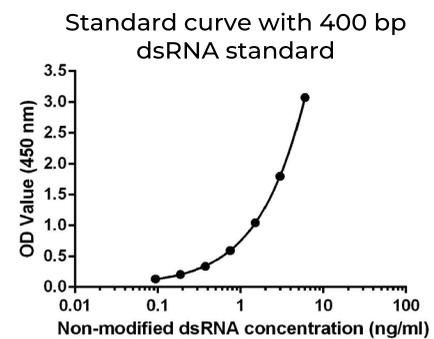


标准曲线



*使用400 bp m1Ψ修饰的dsRNA标准品评估L01020

L01020的线性范围为0.123-30 ng/ml, 灵敏度为0.123 ng/mL,
R2 > 0.99

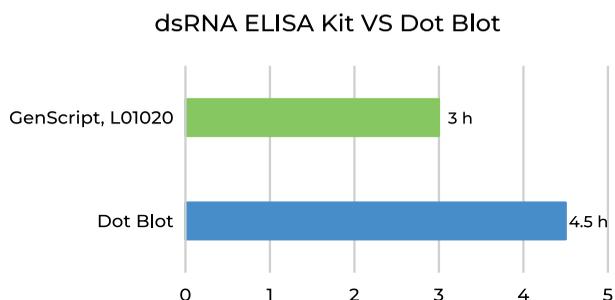


*使用400 bp的非修饰dsRNA标准品评估L01020

L01020的线性范围为0.094-6 ng/ml, 灵敏度为0.094 ng/mL,
R2 > 0.99

产品性能

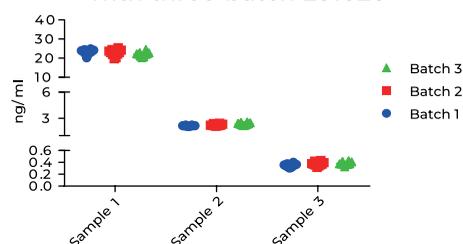
检测时间短



✓ L01020的检测时长为3 h

精密度高

Inter-assay and Intra-assay CV calculation with three-batch L01020



*用3个批次的L01020对3个不同浓度的样本各检测10次，计算变异系数(CV)

✓ L01020的批间批内精密度的CV≤10%

稀释线性佳

Dilution Factor	Expected dsRNA Conc. (ng/mL)	Measured dsRNA Conc. (ng/mL)	Accuracy%
1000	0.738	0.755	102.3
100	0.738	0.787	106.6
50	0.738	0.799	108.3

*使用L01020测试同一样品的不同稀释倍数来评估稀释线性

✓ L01020检测结果不受样本稀释倍数的影响

特异性高

dsRNA Conc. (ng/mL)	mRNA Conc. (ng/mL)	Measured dsRNA Conc. (ng/mL)	Accuracy%
30	300	27.79	92.6
30	30	28.21	94.0
0.123	300	0.138	112.2
0.123	30	0.115	93.5

*两种浓度 (ULOQ 30 ng/mL 和 LLOQ 0.123 ng/mL) 的样品中添加不同量的 mRNA (30 ng/mL 和 300 ng/mL)，以评估 L01020 的特异性

✓ L01020检测结果不受基质的影响



扫码关注公众号

地址: 江苏省南京市江宁科学园雍熙路28号
 邮箱: product@genscript.com.cn
 售后电话: 400-025-8686-6707

DREK20241202

网址: www.genscript.com.cn
 电话: 400-0258686 转 5810
 售后邮箱: cpbu.techsupport@genscript.com

