

Batch information

Tag-lite ligand 2 HTRF D2 receptor red antag. - 5K pts
Product reference : **L0002RED**

Product characteristics :

- Lot Number	07A
- Conditioning buffer	HEPES 50 mM pH 7.4 + 10% DMSO
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2023.12.01

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Pro)	0.36 nmoles	1 295

PLEASE NOTE

The expiration date of batch 07A has been extended to 2023-12-01 after validation by QC.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

