

Batch information

Tag-lite ligand 2

Dopamine D2 receptor red antago. - 5,000 tests

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	2022.05.01

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 2022-05-01 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

