

## Batch information

### Tag-lite ligand 2

HTRF D2 receptor red antag. - 5K pts

Product reference :

**L0002RED**

### Product characteristics :

- Lot Number	07B
- Conditioning buffer	HEPES 50 mM pH 7.4 + 10% DMSO
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2024.03.01

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Pro)	0.49 nmoles	1 295

### PLEASE NOTE

*Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)*

