

Batch information

STD

HTRF CTLA4/CD80 Bind. kit - Std

Product reference :

64CTLA80CDA

Product characteristics :

- Lot Number	02A
- Conditioning buffer	tampon hepes 50mM pH7 + 01% BSA + stabilizer
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2024.11.18

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	0.13 nmoles	N/A

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

