

Batch information

STD

HTRF FLAG check kit - Std

Product reference :

63ADK036CDA

Product characteristics :

- Lot Number	17A
- Conditioning buffer	PO4 50 Mm pH7, BSA 0.1% + stabilizer
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2024.11.23

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ag)	0.50 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

