

Batch information

STD

HTRF Human ApoE Kit - Std

Product reference :

64APOECDA

Product characteristics :

- Lot Number	01A
- Conditioning buffer	TP PO4 50 mM pH 7 + 0.2% BSA + stabilizer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.10.12
- Volume per vial	50.00 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Apo)	3 000.00 ng	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

