

Batch information

phospho AKT-3

HTRF AKT3 p-S473 kit - Ctrl lysate

Product reference :

63ADK082TDA

Product characteristics :

- Lot Number	09A
- Conditioning buffer	LB1
- Appearance	Frozen form
- Storage conditions	-80 °C
- Expiration date	2024.01.18

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	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

