

Batch information

control lysate

HTRF MLKL p-S358 Kit - Ctrl lysate

Product reference :

64MLKLS8TDA

Product characteristics :

- Lot Number	01A
- Conditioning buffer	Lysis buffer
- Appearance	Frozen form
- Storage conditions	-80 °C
- Expiration date	2024.07.12

Weight indications :

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

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

