

Batch information

Control lysate

HTRF EIF2 alpha p-S52 kit - Ctrl lysate

Product reference :

64EF2TDA

Product characteristics :

- Lot Number	05A
- Conditioning buffer	Lysis Buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.12.15
- Volume per vial	150.00 µL

Weight indications :

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

PLEASE NOTE

