

Batch information

Control lysate

VEGFR2 p-Y1175 kit - control lysate

Product reference :

63ADK041TDA

Product characteristics :

- Lot Number	12A
- Conditioning buffer	Lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2023.12.03

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	1

PLEASE NOTE

