

Batch information

control lysate

HTRF FGFR1 p-Y653/654 kit - Ctrl lysate

Product reference :

64FGFR1Y6TDA

Product characteristics :

- Lot Number	03A
- Conditioning buffer	Lysis buffer
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2024.11.24

Weight indications :

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

PLEASE NOTE

