

Batch information

phosph MEK1/2-K

CUSTOM LABELING

Product reference :

61P17KAY

Product characteristics :

- Lot Number	31A
- Packaging Unit	100 000 tests/vial
- Conditioning buffer	Hépès 0.05M pH7 + 0.1%BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2031.10.06
- Volume per vial	2.50 mL
- Active Moiety Concentration (Ab)	120.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	300.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	3.00 ng
Number of tests/vial	100 000

*Actual amount per well should be optimized according to assay conditions

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

