

Tracer Dasatinib-Red - 20 nmol
Product reference : **62KB02REDE**

Product characteristics :

- Lot Number _____ 02A
- Conditioning buffer _____ -
- Appearance _____ Frozen form
- Storage conditions _____ <-16 °C
- Expiration date _____ 2023.02.19

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Kin)	800.00 µl	N/A

PLEASE NOTE

The expiration date of batch 02A has been extended to 2023-02-19 after validation

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

