

D2V2-NHS

d2v2-NHS - 0.1 mg

Product reference :

65DVSABA

Product characteristics :

- Lot Number_____ 09A
- Conditioning buffer_____ ///
- Appearance_____ Desiccated form
- Storage conditions_____ <-16 °C
- Expiration date_____ 2024.10.19

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (DV)	0.10 mg	1 066

PLEASE NOTE

Rz (D0650nm/280nm)=34.7

Reconstitute with DMSO

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