

Product Information

Description:	Activated DNA for PARP assay
Applications:	Study enzyme kinetics, and screen small molecular inhibitors of PARP for drug discovery and HTS applications.
Supplied As:	Aqueous buffer solution
Quantity:	500 µl
Storage:	-20°C or -80°C