

Product Information

Description:	PARP Substrate Mixture 2 is a mix of biotinylated and non-biotinylated substrate (NAD ⁺).
Supplied As:	Tris-buffered solution.
Storage:	Upon receipt, store at -80°C.
Stability:	When stored at -80°C, the product is stable for six months after first thaw. Avoid repeated freeze/thaw cycles.
Application:	<p>It is suitable for use in PARP assays using cell extracts or purified proteins. This mixture has been optimized for use:</p> <ul style="list-style-type: none">• PARP 6 Chemiluminescent Assay Kit (BPS Bioscience #80556)• PARP 7 Chemiluminescent Assay Kit (BPS Bioscience #79729)• PARP 10 Chemiluminescent Assay Kit (BPS Bioscience #80560)• PARP 11 Chemiluminescent Assay Kit (BPS Bioscience #80561)• PARP 14 Chemiluminescent Assay Kit (BPS Bioscience #80568)• PARP 15 Chemiluminescent Assay Kit (BPS Bioscience #80567)• TNKS1 (PARP5A) Chemiluminescent Assay Kit (BPS Bioscience #78405)• TNKS2 (PARP5B) Chemiluminescent Assay Kit (BPS Bioscience #78406) <p>Use in other assays may require optimization.</p>