

Product Information

Description:	PARP Substrate Mixture 1 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 1 Chemiluminescent Assay Kit (BPS Bioscience #80551)• PARP 2 Chemiluminescent Assay Kit (BPS Bioscience #80552)• PARP 3 Chemiluminescent Assay Kit (BPS Bioscience #80553) <p>Use in other assays may require optimization.</p>