

## Product Information

**Description**

A set of Poly (ADP-ribose) polymerases (PARP) inhibitors (PARPi) including PARPi with a broad binding specificity toward PARPs like Olaparib, Niraparib, Rucaparib, Talazoparib, Veliparib, as well as more selective PARPi like AZD5305 specific toward PARP1, XAV939 specific toward TNSK1 and TNSK2 and RBN-2397 specific toward PARP7.

**Concentration:**

10 mM solution

**Format:**

DMSO solution

**Stability:**

At least 6 months at -80°C. Avoid freeze/thaw cycles.

**Storage:**

-80°C

**Optimized for Use With:**

- PARP Chemiluminescent Assay Kits (BPS Bioscience #80569, 80551, 80552, 80553, 80573, 80579, 80578, 80572, 80556, 79729, 80560, 80561, 80568 and 80567)
- PARPtrap™ Assay Kits for PARP1 and for PARP2 (BPS Bioscience #80584 and #78296)
- PARPtrap™ Combo Assay Kit for PARP1 and PARP2 (BPS Bioscience #78317)