

Product Information

Description:	The ADP-Ribosylation Cycle Inhibitor Mix is a 500x cocktail of two inhibitors (AZD2281 and PDD00017273). AZD2281 (Olaparib) is a PARP inhibitor with broad pan-PARP activity, which blocks a large proportion of the ADP-ribosylation of proteins. PDD00017273 is a potent, selective poly (ADP-ribose) glycohydrolase (PARG) inhibitor that blocks PARG-mediated de-ribosylation of proteins. Used together, these two inhibitors “freeze” the state of cellular ribosylation at the time the mix is applied to the cells. This product is designed to supplement a cell lysis buffer of choice to prevent post-lysis changes in the ADP-ribosylation status of extracted proteins.
Formulation:	5 mM AZD2281 (Olaparib) (BPS Bioscience #27003) and 5 mM PDD00017273 in 100% DMSO (Dimethyl sulfoxide).
CAS Registry:	Olaparib: 763113-22-0 PDD00017273: 1945950-21-9
Formula:	Olaparib: $C_{24}H_{23}FN_4O_3$ PDD00017273: $C_{23}H_{26}N_6O_4S$
MW:	Olaparib: 434.46 PDD00017273: 514.62
Storage/Stability:	Upon receipt store at $-80^{\circ}C$. Stable for 6 months from date of receipt, when stored as directed.
Instructions for Use:	Thaw right before use. Dilute 500-fold directly in the cell lysis buffer of choice. Aliquots are single use and should be discarded after thawing. Discard excess lysis buffer containing the inhibitor mix at the end of the day. <i>Note: This inhibitor mix is designed for the analysis of proteins that are ribosylated by PARP1/2/3/4 and PARP10 (approximately 80% of proteins) and are de-ribosylated by PARG. For proteins that are the target of other PARPs or a ribosylation eraser other than PARG, different inhibitors may be necessary.</i>
Shipping Temperature:	$-80^{\circ}C$.