

PARP Screening Services

Poly (ADP-ribose) Polymerase (PARP) plays important roles in chromatin structure, DNA repair, and apoptosis. PARP is overexpressed in a number of human cancers, and its expression correlates with an aggressive phenotype and resistance to DNA damage-inducing therapeutic agents. PARP inhibitors are an exciting new class of anti-cancer drugs that show particular promise for breast and ovarian cancers with mutations in the *PTEN*, *BRCA1* or *BRCA2* genes.

There are at least 16 different members of the PARP family, which vary widely in their function, tissue distribution, and subcellular localization. BPS Bioscience offers a unique panel of 12 different PARP isozymes, including tankyrase 1 (PARP5A) and tankyrase 2 (PARP5B) for evaluation of lead compounds.

PARP Assays		
PARP1	PARP5B (TNKS2)	PARP10
PARP2	PARP6	PARP11
PARP3	PARP7	PARP12
PARP5A (TNKS1)	PARP8	PARP14
		PARP15

*To review sample data for the PARP assays listed above, click on the following link or copy and paste it into your web browser: http://www.bpsbioscience.com/images/pdf/PARP_profiling_data.pdf

- Fast turn-around time, typically under 2 weeks
- Determine IC50 values (10 point, duplicate) or screen a single concentration over our full panel of isozymes.
- Standard enzyme panels available, or customize to meet your needs
- Enzyme activity is validated on every assay to ensure reliable results
- Choice of dose range and concentrations

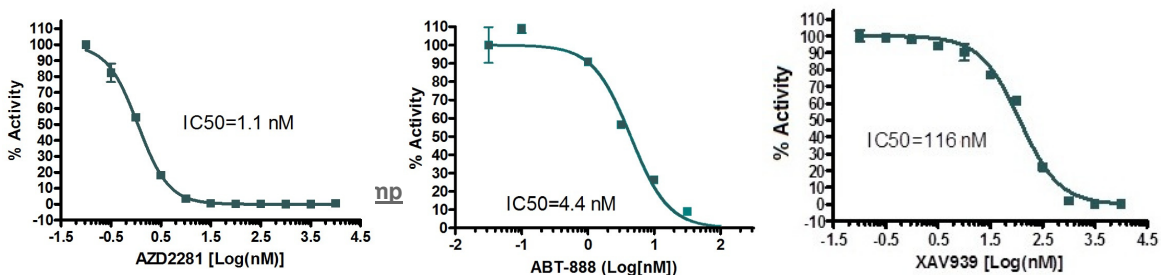


Figure 1. Inhibition of PARP1 activity by AZD2281, ABT-888 and XAV-939 using BPS Bioscience's PARP screening and profiling services.

Contact us for a free quotation or information on our PARP Screening and Profiling Services by visiting:

www.bpsbioscience.com/contact-us