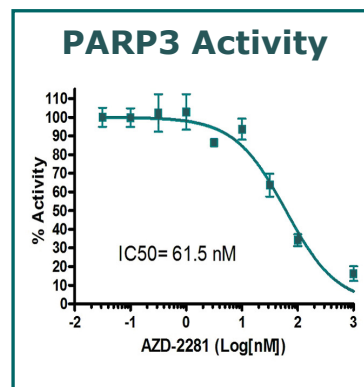
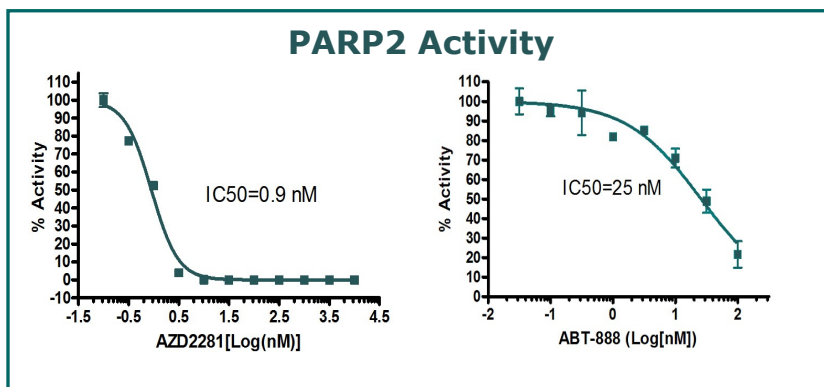
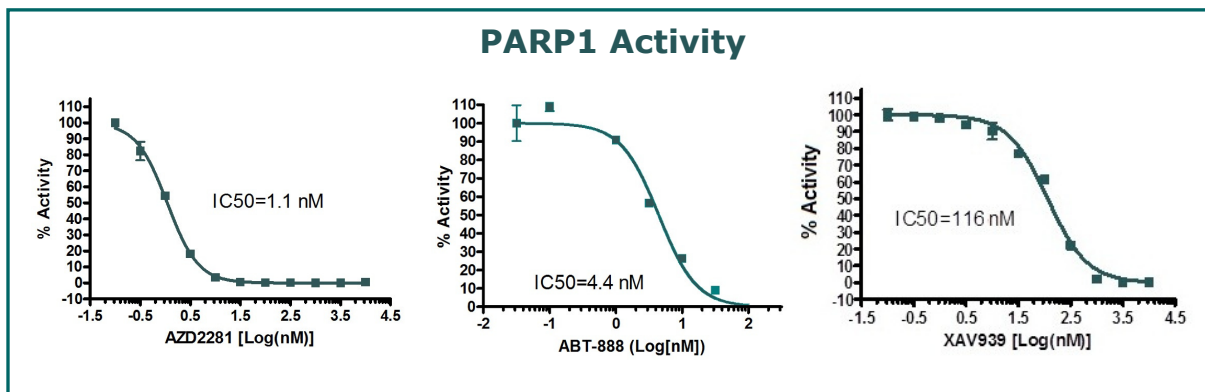


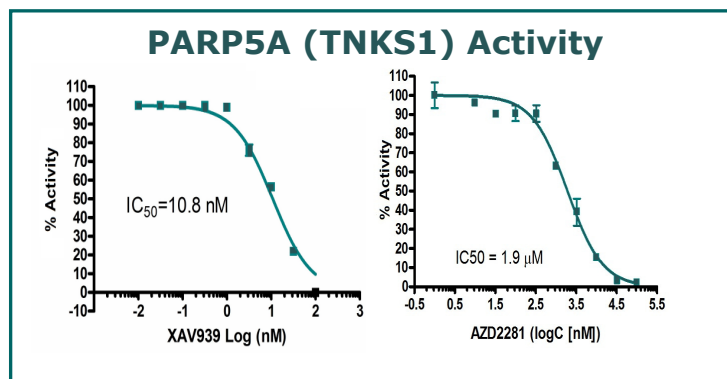
Poly (ADP-ribose) Polymerase (PARP) Profiling Services



Only from BPS Bioscience!

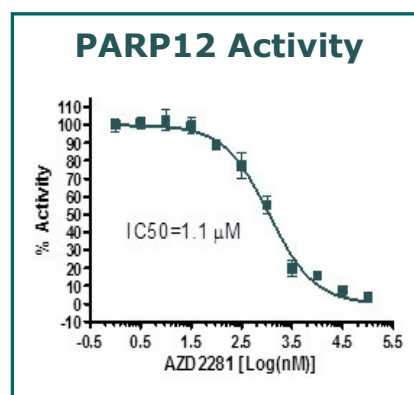
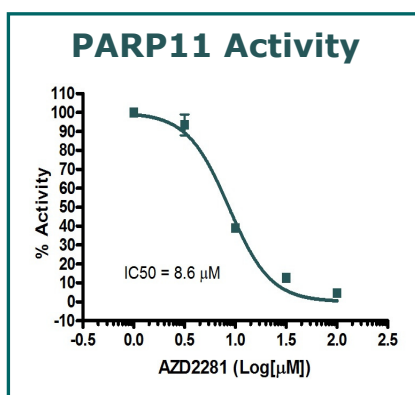
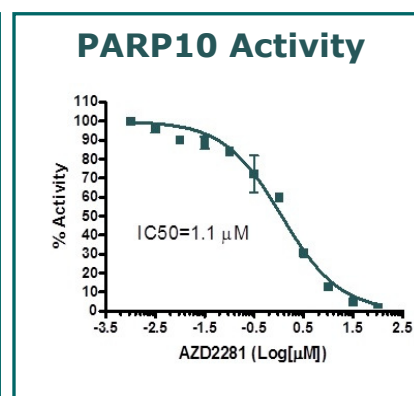
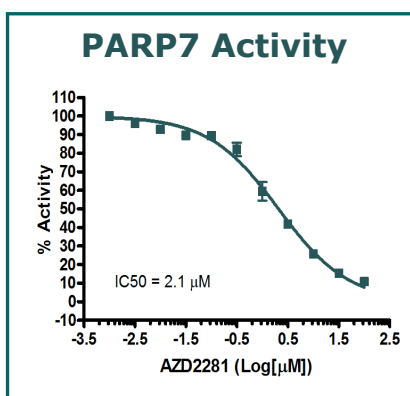
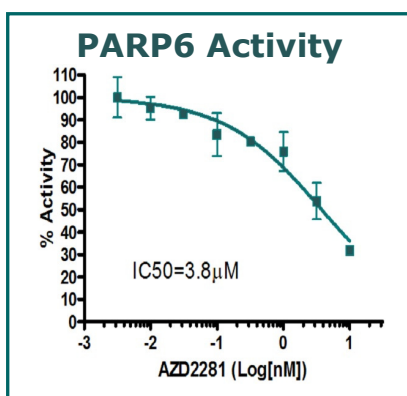
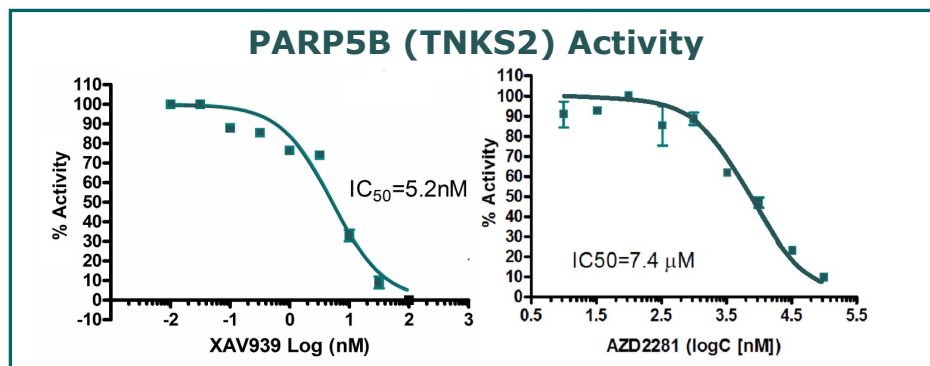
Profiling and Screening of PARP inhibitors or activators. Our panel includes 10 different PARP isozymes: PARPs 1, 2, 3, 5A (Tankyrase I), 5B (Tankyrase II), 6, 7, 10, 11, and 12.

Contact us for a quote today!



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
 To place your order, please contact us by Phone **1.858.829.3082**, Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



Also available: IC₅₀s for **HDACs 1-11**, **DNMTs (1, 3A, & 3B)**, **demethylases (LSD1, JMJD2A-C)**, and 15 different histone **methyltransferases**, including **EZH2**, **MLL**, **G9a**, and **PRMT4 (CARM1)**.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082**, Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com