

Bromodomain Screening Services

Acetylation of lysine residues is a dynamic process that directs gene transcription and chromatin structure. Bromodomain (BRD)-containing proteins recognize and bind to these acetylated lysines to regulate epigenetically controlled processes. Recent research has implicated BRDs in a number of human diseases, including diabetes, cancer, and inflammatory disorders. The identification and optimization of small molecule inhibitors is essential as BRDs emerge as an important new drug target.

BPS Bioscience offers a broad panel of 25 different Bromodomains for evaluation of lead compounds.

Bromodomain Assays			
ATAD2A	BRD2 (BD1 + BD2)	BRD4 (BD1 + BD2)	CREBBP
ATAD2B	BRD3 (BD1)	BRD9 (BD1)	SMARCA2
BAZ2B	BRD3 (BD2)	BRDT (BD1)	TAF1 (BD2)
BRD1	BRD3 (BD1 + BD2)	BRG1	TAF1 (BD1 + BD2)
BRD2 (BD1)	BRD4 (BD1)	BRPF3	TAF1L (BD2)
BRD2 (BD2)	BRD4 (BD2)	CECR2	TAF1L (BD1 + BD2)
			TRIM24 (TIF1)

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

Sample Results: JQ1 Inhibition Curves

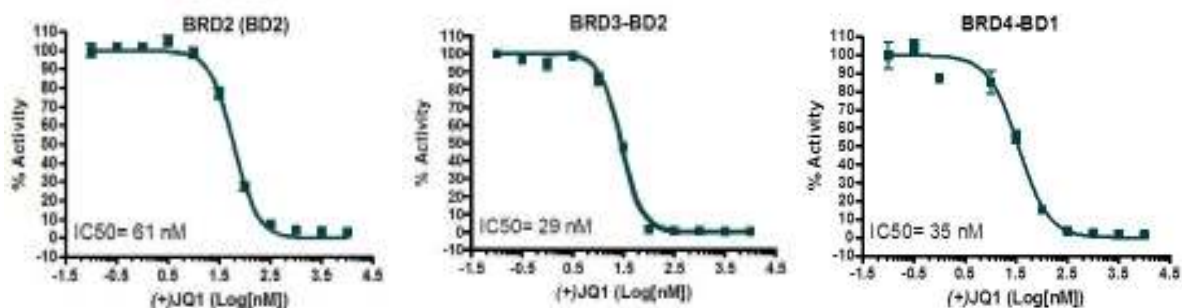


Figure 1. Inhibition of BRD2-BD2, BRD3-BD2 and BRD4-BD1 by (+)-JQ1 using BPS Bioscience's Bromodomain screening and profiling services, proteins and inhibitors.

Contact us for a free quotation or information on our Bromodomain Screening and Profiling Services by visiting: www.bpsbioscience.com/contact-us