



## Bromodomains

Bromodomains (BRDs) are protein domains which recognize acetylated lysines. Bromodomains are emerging as a new class of targets, alongside HATs and HDACs, which will be important in cancer drug discovery as well as treatments of inflammatory diseases such as cardiovascular disease and rheumatoid arthritis. There are over 50 proteins containing one or more BRDs. BRDs have been identified in HATs, methyltransferases (e.g. MLL), chromatin-remodeling complexes, and transcriptional coactivators.

BPS Bioscience provides an ever-growing portfolio of BRD enzymes, assay kits, and inhibitors suitable for bromodomain profiling, inhibitor screening assays, and drug discovery.

### BRD Inhibitors

Inhibitor	Cat. #	Size
(+) - JQ1	27400	10 mg
	27401	1 mg
	27402	5 mg
RediSolution™ (+)-JQ1	27403	1 mg
(-)-JQ1	27334	5 mg
	27335	25 mg
CHS-828	27333	25 mg
PFI-1	27332	25 mg
RVX-208	27111	5 mg
I-CBP112 (Hydrochloride)	27336	1 mg
I-BET151 (GSK1210151A)	27629-1	5 mg
	27629-2	10 mg
	27629-3	50 mg
I BET-762	27630-1	5 mg
	27630-2	10 mg
OTX-015	27636-1	5 mg
	27636-2	10 mg
	27636-3	25 mg
SGC-CBP30	27650-1	10 mg
	27650-2	50 mg

### BRD Screening/Profiling Kits

AlphaScreen® Assay Kit	Cat. #	Size
ATAD2A Inhibitor Screening Assay Kit	32601	384 rxns
ATAD2B Inhibitor Screening Assay Kit	32605	384 rxns
BAZ2B Inhibitor Screening Assay Kit	32600	384 rxns
BRD1 Inhibitor Screening Assay Kit	32521	384 rxns
BRD2 (BD2) Inhibitor Screening Assay Kit	32522	384 rxns
BRD3 (BD1) Inhibitor Screening Assay Kit	32513	384 rxns
BRD3 (BD2) Inhibitor Screening Assay Kit	32523	384 rxns
BRD4 (BD1) Inhibitor Screening Assay Kit	32514	384 rxns
BRD4 (BD2) Inhibitor Screening Assay Kit	32524	384 rxns
BRD9 (BD1) Inhibitor Screening Assay Kit	32519	384 rxns
SMARCA4/BRG1 Inhibitor Screening Assay Kit	32609	384 rxns
BRPF3 Inhibitor Screening Assay Kit	32608	384 rxns
CREBBP Inhibitor Screening Assay Kit	32607	384 rxns
SMARCA2 Inhibitor Screening Assay Kit	32610	384 rxns
TAF1 (BD1+BD2) Inhibitor Screening Assay Kit	32604	384 rxns
TAF1 (BD2) Inhibitor Screening Assay Kit	32624	384 rxns
TAF1L (BD1+BD2) Inhibitor Screening Assay Kit	32603	384 rxns
TAF1L (BD2) Inhibitor Screening Assay Kit	32602	384 rxns
TRIM24 Inhibitor Screening Assay Kit	32606	384 rxns

TR-FRET Assay Kit	Cat. #	Size
ATAD2A TR-FRET Assay Kit	32618	384 rxns
ATAD2B TR-FRET Assay Kit	32620	384 rxns
BRD1 TR-FRET Assay Kit	32625	384 rxns
BRD2 (BD1+BD2) TR-FRET Assay Kit	32614	384 rxns
BRD3 (BD1+BD2) TR-FRET Assay Kit	32615	384 rxns
BRD4 (BD1) TR-FRET Assay Kit	32613	384 rxns
BRD4 (BD2) TR-FRET Assay Kit	32617	384 rxns
BRD4 (BD1+BD2) TR-FRET Assay Kit	32612	384 rxns
BRDT (BD1+BD2) TR-FRET Assay Kit	32616	384 rxns
BRD9 TR-FRET Assay Kit	32621	384 rxns
BRPF3 TR-FRET Assay Kit	32626	384 rxns
BPTF/FALZ TR-FRET Assay Kit	32632	384 rxns
CECR2 TR-FRET Assay Kit	32622	384 rxns
CREBBP TR-FRET Assay Kit	32619	384 rxns
TAF1 (BD2) TR-FRET Assay Kit	32627	384 rxns
TAF1 (BD1+BD2) TR-FRET Assay Kit	32628	384 rxns
TAF1L (BD1+BD2) TR-FRET Assay Kit	32629	384 rxns
TRIM24 TR-FRET Assay Kit	32630	384 rxns
WDR9/BRWD1 TR-FRET Assay Kit	32631	384 rxns

## Bromodomain Proteins

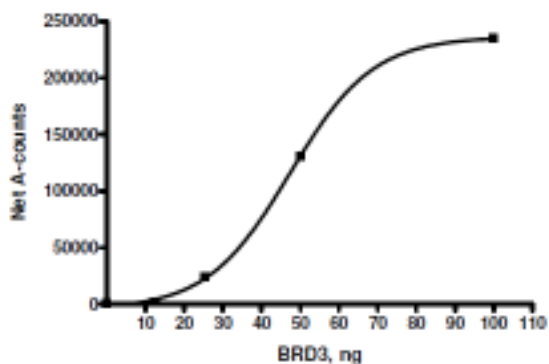
Protein	Cat. #	Size
ASH1L, His-tag	31147	100 µg
ATAD2A, His-tag	31109	100 µg
ATAD2A, GST-tag	31121	100 µg
ATAD2B, His-tag	31117	100 µg
ATAD2B, GST-tag	31125	100 µg
BAZ1A, GST-tag	31136	100 µg
BAZ1B, GST-tag	31137	100 µg
BAZ1B, His-tag	31145	100 µg
BAZ2B, His-tag	31113	100 µg
BRD, His-tag	31010	100 µg
BRD1, GST-tag	31011	100 µg
BRD2 (BD1), GST-tag	31021	100 µg
BRD2 (BD1), His-tag	31022	100 µg
BRD2 (BD2), GST-tag	31023	100 µg
BRD2 (BD2), His-tag	31020	100 µg
BRD2 (BD1+BD2), GST-tag	31024	100 µg
BRD2 (BD1+BD2), His-tag	31025	100 µg
BRD3 (BD1), GST-tag	31032	100 µg
BRD3 (BD1), His-tag	31030	100 µg
BRD3 (BD2), GST-tag	31033	100 µg
BRD3 (BD2), His-tag	31031	100 µg
BRD3 (BD1+BD2), His-tag	31034	100 µg
BRD3 (BD1+BD2), GST-tag	31035	100 µg
BRD4 (BD1), GST-tag	31040	100 µg
BRD4 (BD1), His-tag	31042	100 µg
BRD4 (BD2), GST-tag	31041	100 µg
BRD4 (BD2), His-tag	31043	100 µg
BRD4 (BD1+BD2), GST-tag	31044	100 µg
BRD4 (BD1+BD2), His-tag	31045	100 µg
BRD4 (BD1+BD2), un-tagged	31047	100 µg
BRD9, His-tag	31090	100 µg
BRD9, GST-tag	31091	100 µg
BRDT (BD1), His-tag	31101	100 µg
BRDT (BD1), GST-tag	31108	100 µg
BRDT (BD2), His-tag	31100	100 µg
BRDT (BD2), GST-tag	31060	100 µg
BRDT (BD1+BD2), GST-tag	31061	100 µg

Protein	Cat. #	Size
BRPF1, His-tag	31112	100 µg
BRPF3, GST-tag	31130	100 µg
CECR2, GST-tag	31138	100 µg
CECR2, His-tag	31046	100 µg
CREBBP, His-tag	31119	100 µg
CREBBP, GST-tag	31128	100 µg
FALZ (BPTF), GST-tag	31134	100 µg
FALZ (BPTF), His-tag	31131	100 µg
GCN5 (727-837), His-tag	31114	100 µg
P300 (1046-1163), His-tag	31118	100 µg
PB1 (BD4), His-tag	31122	100 µg
PB1 (BD6), His-tag	31133	100 µg
PCAF (KAT2B), His-tag	31120	100 µg
PHIP (BD1), GST-tag	31142	100 µg
PHIP (BD1), His-tag	31140	100 µg
SMARCA2, GST-tag	31141	100 µg
SMARCA2, His-tag	31111	100 µg
SMARCA4/BRG1 (BD1), GST-tag	31132	100 µg
SMARCA4/BRG1 (BD1), His-tag	31102	100 µg
TAF1 (BD1), His-tag	31123	100 µg
TAF1 (BD2), GST-tag	31126	100 µg
TAF1 (BD2), His-tag	31110	100 µg
TAF1 (BD1+BD2), GST-tag	31124	100 µg
TAF1L (BD1), GST-tag	31105	100 µg
TAF1L (BD1), His-tag	31103	100 µg
TAF1L (BD1+BD2), GST-tag	31107	100 µg
TAF1L (BD2), His-tag	31104	100 µg
TAF1L (BD2), GST-tag	31106	100 µg
TRIM24 (TIF1), His-tag	31116	100 µg
TRIM24 (TIF1), GST-tag	31127	100 µg
TRIM28, GST-tag	31139	100 µg
TRIM28, His-tag	31146	100 µg
WDR9/BRWD1 (BD2), GST-tag	31135	100 µg
WDR9/BRWD1 (BD2), His-tag	31115	100 µg
ZMYND8, GST-tag	31143	100 µg
ZMYND11, GST-tag	31144	100 µg

## BRD-Related Products

Product	Cat. #	Size
3x ATAD2 Detection Buffer	33006	10 ml
3x ATAD2 Assay Buffer	33007	10 ml
3x BRD assay buffer	33001	10 ml
3x BRD detection buffer	33002	10 ml
BET Bromodomain Ligand	33000	500 µl
Bromodomain Ligand 2	33003	500 µl
Bromodomain Ligand 3	33008	500 µl
Bromodomain Ligand 4	33010	500 µl
Bromodomain Non-acetylated Ligand 1	33005	500 µl
Bromodomain Non-acetylated Ligand 2	33004	500 µl
Bromodomain Non-acetylated Ligand 3	33009	500 µl
Bromodomain Non-acetylated Ligand 4	33011	500 µl

### BRD3 (BD1) binding to the acetylated substrate



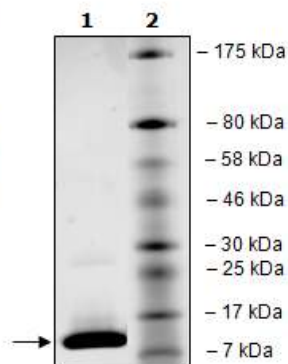
BRD3 (BD1) binding activity of BRD3 (29-145), BD1 (Cat. #31030) measured using the BRD3 (BD1) Inhibitor Screening Assay Kit, Cat. #32513.

### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
6.4 µg BRD4

**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 14.7 kDa  
**Purity:** ≥98%



Purity and gel image of BRD4 (342 – 460), Cat. # 31043.

6044 Cornerstone Ct. West  
Suite E  
San Diego, CA 92121  
Ph: 858.829.3082  
Fax: 858.481.8694  
[www.bpsbioscience.com](http://www.bpsbioscience.com)