

## SPR study of BRD4-1

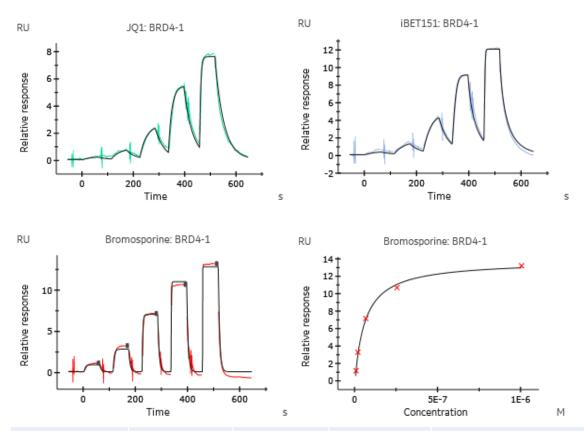
(Bromodomain-containing protein 4 domain 1, Protein HUNK1)

CAT#: BRD4-1

Enzyme: Human BRD4-1

Scientific Information: O60885 (<a href="http://www.expasy.org/uniprot/O60885">http://www.expasy.org/uniprot/O60885</a>)

• Close affinity, different kinetics: JQ1 and Bromsporine both showed sub-nano molar binding affinity to the BRD4-1. Bromosporine dissociates about10 times faster than JQ1, which reflects a short on-target residence time.

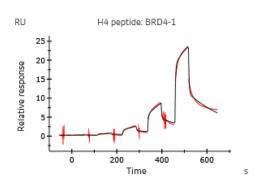


				KD (M) by steady state affinity fitting	to 50% dissociation
Compound	ka (1/Ms)	kd (1/s)	KD (M)		(t <sub>1/2</sub> )
JQ1	1.56E+06	2.14E-02	1.37E-08	N/A	32.4s
Bromosporine	3.15E+06	1.85E-01	5.86E-08	6.45E-8	3.8s



## Reveal Bivalent binder

Data from H4 peptide (Histone H4 peptide (1-21) K5/8/12/16Ac) could only be fitted well to a bivalent analyte model. This agrees with previously published structural studies. As KD1 is significantly lower than KD2, it can be seen as the apparent KD for comparing with IC50 data.



Compound	ka1 (1/Ms)	kd1 (1/s)	ka2 (1/RUs)	kd2 (1/s)	KD1 (M)	KD2 (M)
H4 peptide	3.94E+04	3.46E-01	1.85E-04	2.53E-03	8.78E-06	1.36E+01

## Data reproducibility

Kinetic fitting parameters of three independent run of JQ1 binding to BRD4-1

Replicate	ka (1/Ms)	kd (1/s)	KD (M)
Run 1	1.56E+06	2.14E-02	1.37E-08
Run 2	1.89E+06	2.91E-02	1.54E-08
Run 3	2.08E+06	2.72E-02	1.33E-08

