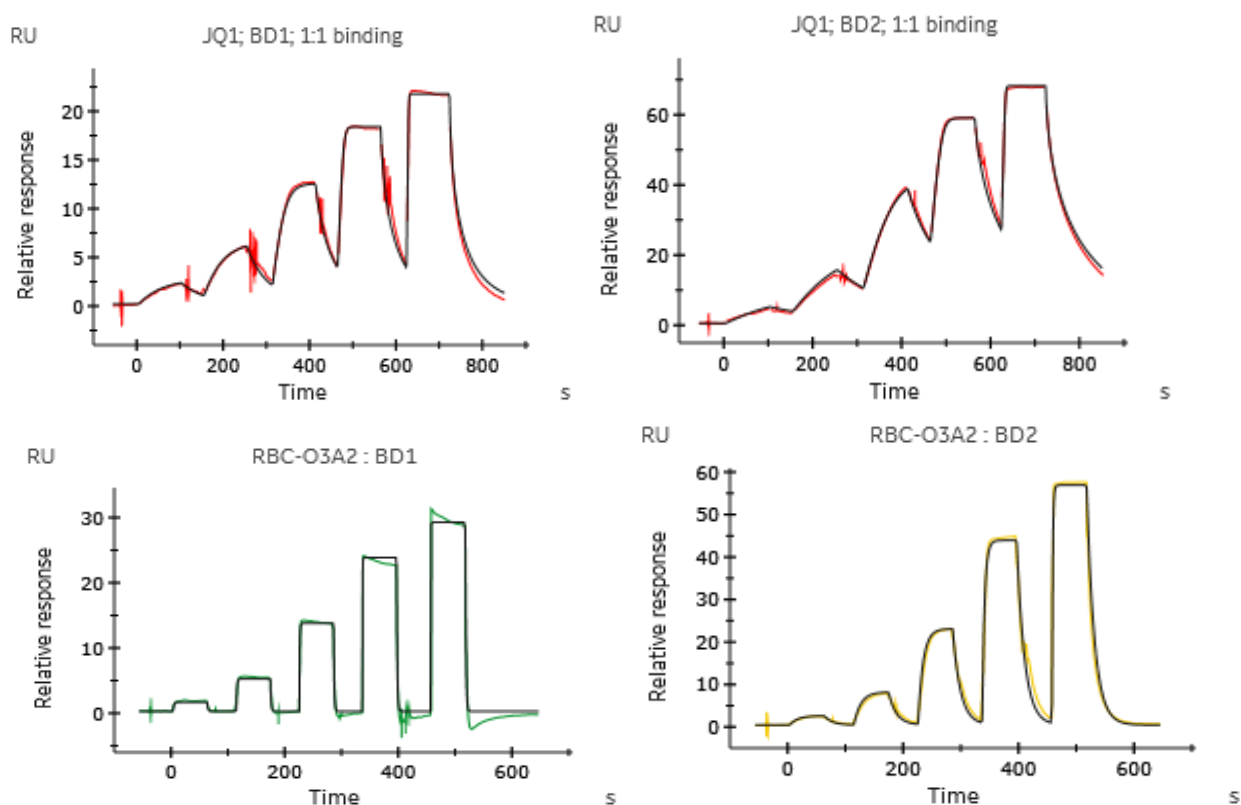


## SPR study of BRD4-2

(Bromodomain-containing protein 4 domain 2, Protein HUNK1)

- **CAT#:** BRD4-2
- **Enzyme:** Human BRD4-2
- **Scientific Information:** O60885 (<http://www.expasy.org/uniprot/O60885>)

**Domain selectivity determined by kinetic profile:** JQ1 and the internal compound RBC-O3A2 were tested against domain 1 (BD1) and domain 2 (BD2) of BRD4 in parallel.



Compound: domain	$k_a$ (1/Ms)	$k_d$ (1/s)	KD (M)	to 50% dissociation ( $t_{1/2}$ ) (s)
JQ1: BD1	1.45E+06	3.04E-02	2.09E-08	22.80
JQ1: BD2	7.82E+05	1.39E-02	1.77E-08	49.86
RBC-O3A2: BD1	5.06E+05	8.37E-01	1.65E-06	0.83
RBC-O3A2: BD2	3.35E+05	7.40E-02	2.21E-07	9.36