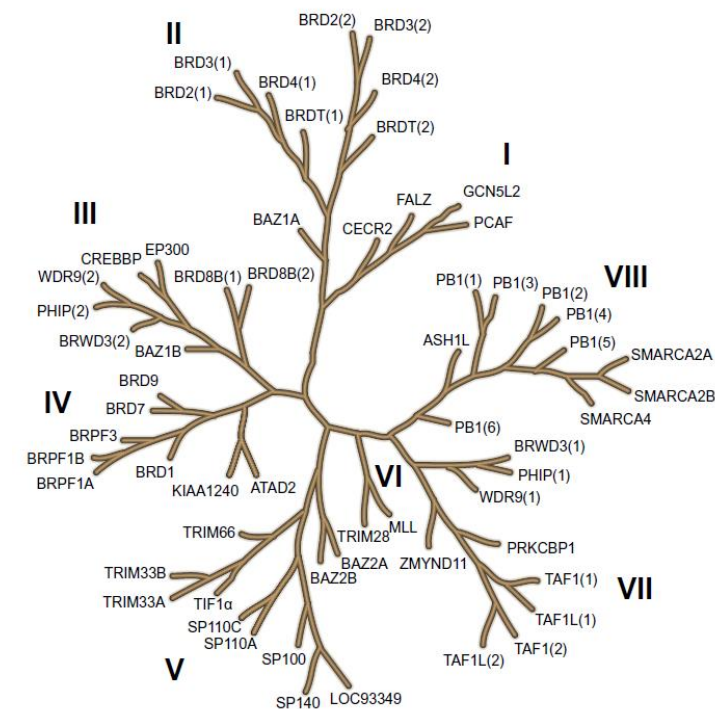




BromoMELTSM

- Extensive coverage of bromodomain family
- Test selectivity of any compound within a few hours
- Available as a service or in kit format
- Fast and easy to perform in your own lab

	1	2	3	4	5	6	7	8	9	10	11	12
A	color reference	Buffer blank	BRD2-1 (His)	BRD2-2 (His)	BRD2-Tndm (His)	ATAD2 (GST)	ATAD2 (His)	ATAD2B (His)	TAF1-1 (His)	TAF1-2 (GST)	TAF1-2 (His)	
B		CECR2 (GST)	BRD3-1 (His)	BRD3-2 (His)	BRD3-Tndm (His)	BRD1 (His)	BRD7 (GST)	BRD9 (His)	TAF1L-1 (GST)	TAF1L-1 (His)	TAF1L-2 (His)	
C		BPTF-[PHD-BRD] (His)	BRD4-1 (His)	BRD4-2 (His)	BRD4-Tndm (His)	BRPF1a (His)	BRPF1b (His)	BRPF3 (His)	TAF1L-Tndm (GST)	TAF1-Tndm (His)	PHIP-Tndm (GST)	
D		BPTF-[PHD-BRD](GST)	BRD4-1 (GST)	BRD4-2 (GST)	BRDT-2 (GST)	BAZ2A (His)	BAZ2B (His)	SP100 (GST)	ASH1L-[BRD](GST)	PB1-1 (GST)	PB1-2 (His)	
E		BPTF-[BRD](His)	BRDT-1 (His)	BRDT-2 (His)	BRDT-Tndm (His)	SP100 (His)	SP110c (GST)	SP140 (His)	PB1-3 (His)	PB1-4 (His)	PB1-5 (His)	
F		KAT2A (His)	BAZ1B (GST)	BAZ1B (His)	BRD8-1 (GST)	SP140L (His)	TRIM24 (His)	Buffer blank	PB1-6 (GST)	SMARCA2a (GST)	SMARCA2a (His)	
G		KAT2B (His)	BRD8-1 (his)	BRWD1-2 (His)	BRWD3-2 (GST)	TRIM33a (GST)	TRIM33a (His)	TRIM28 (GST)	SMARCA2 b (GST)	SMARCA4 (GST)	SMARCA4 (His)	
H		KAT2B (GST)	CREBBP (His)	EP300 (His)	PHIP-2 (His)	TRIM33b (His)	TRIM66 (His)	TRIM28 (His)	BAZ1A (His)	Buffer	GST alone	
		Subfamily 1	Subfamily 2	Subfamily 3		Subfamily 4	Subfamily 5	Subfamily 6	Subfamily 7	Subfamily 8		



Filippakopoulos et al, FEBS Letters 586 (2012) 2692-2704



BromoMELTSM

Sample data

Subfamily 2

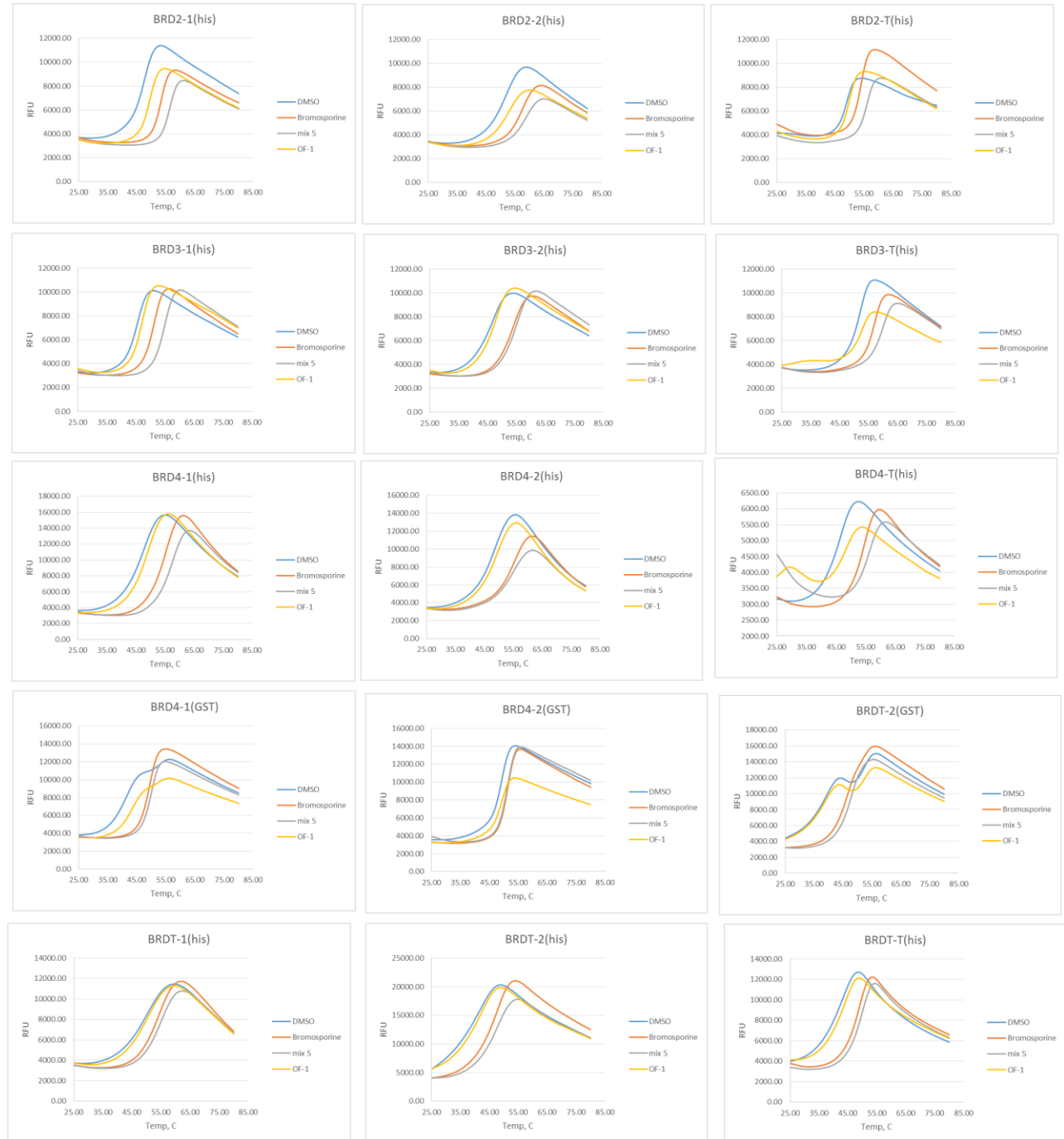
BRD2-1 (His)	BRD2-2 (His)	BRD2-Tndm (His)
BRD3-1 (His)	BRD3-2 (His)	BRD3-Tndm (His)
BRD4-1 (His)	BRD4-2 (His)	BRD4-Tndm (His)
BRD4-1 (GST)	BRD4-2 (GST)	BRDT-2 (GST)
BRDT-1 (His)	BRDT-2 (His)	BRDT-Tndm (His)

Subfamily 2

		Tm	dTm
BRD2-1 (His)	DMSO	47.5	0
BRD2-1 (His)	Bromosporine	53.5	6
BRD2-1 (His)	mix 5	56.5	9
BRD2-1 (His)	OF-1	49.5	2
BRD2-2 (His)	DMSO	52	0
BRD2-2 (His)	Bromosporine	57.5	5.5
BRD2-2 (His)	mix 5	59	7
BRD2-2 (His)	OF-1	53	1
BRD2-Tndm (His)	DMSO	49	0
BRD2-Tndm (His)	Bromosporine	54	5
BRD2-Tndm (His)	mix 5	56	7
BRD2-Tndm (His)	OF-1	50	1
BRD3-1 (His)	DMSO	45.5	0
BRD3-1 (His)	Bromosporine	51.5	6
BRD3-1 (His)	mix 5	55	9.5
BRD3-1 (His)	OF-1	47	1.5
BRD3-2 (His)	DMSO	47	0
BRD3-2 (His)	Bromosporine	54	7
BRD3-2 (His)	mix 5	55.5	8.5
BRD3-2 (His)	OF-1	48	1
BRD3-Tndm (His)	DMSO	52.5	0
BRD3-Tndm (His)	Bromosporine	57	4.5
BRD3-Tndm (His)	mix 5	60	7.5
BRD3-Tndm (His)	OF-1	53	0.5
BRD4-1 (His)	DMSO	48	0
BRD4-1 (His)	Bromosporine	55.5	7.5
BRD4-1 (His)	mix 5	57.5	9.5
BRD4-1 (His)	OF-1	49	1
BRD4-2 (His)	DMSO	48.5	0
BRD4-2 (His)	Bromosporine	55.5	7
BRD4-2 (His)	mix 5	55	6.5
BRD4-2 (His)	OF-1	49.5	1

		Tm	dTm
BRD4-Tndm (His)	DMSO	46.5	0
BRD4-Tndm (His)	Bromosporine	54.5	8
BRD4-Tndm (His)	mix 5	56.5	10
BRD4-Tndm (His)	OF-1	48.5	2
BRD4-1 (GST)	DMSO	41.5	0
BRD4-1 (GST)	Bromosporine	49.5	8
BRD4-1 (GST)	mix 5	50	8.5
BRD4-1 (GST)	OF-1	43.5	2
BRD4-2 (GST)	DMSO	49.5	0
BRD4-2 (GST)	Bromosporine	51.5	2
BRD4-2 (GST)	mix 5	51.5	2
BRD4-2 (GST)	OF-1	49.5	0
BRDT-1 (His)	DMSO	50	0
BRDT-1 (His)	Bromosporine	54.5	4.5
BRDT-1 (His)	mix 5	55.5	5.5
BRDT-1 (His)	OF-1	50.5	0.5
BRDT-2 (His)	DMSO	42	0
BRDT-2 (His)	Bromosporine	48	6
BRDT-2 (His)	mix 5	48	6
BRDT-2 (His)	OF-1	42.5	0.5
BRDT-2 (GST)	DMSO	39.5	0
BRDT-2 (GST)	Bromosporine	46.5	7
BRDT-2 (GST)	mix 5	48	8.5
BRDT-2 (GST)	OF-1	39	-0.5
BRDT-Tndm (His)	DMSO	43.5	0
BRDT-Tndm (His)	Bromosporine	49	5.5
BRDT-Tndm (His)	mix 5	50	6.5
BRDT-Tndm (His)	OF-1	44	0.5

mix 5: JQ1/CBP112/IBET151/GKS2801/PFI3
Compound concentration in assay: 25 μ M





BromoMELTSM

Sample data: compound fingerprints

