

# PRODUCT INFORMATION



## Bromodomain Inhibitor Set Item No. 28315

Page 1 of 2

The Bromodomain Inhibitor Set consists of approximately 35 selective and non-selective bromodomain inhibitors and negative controls. This set contains bromodomain inhibitors targeting members of the bromodomain and extra-terminal domain (BET) and bromodomain and plant homeodomain finger-containing (BRPF) families, CBP/CREBBP and p300, BRD9 and BRD7, and TAF1, among others. The composition of this inhibitor set may vary somewhat depending upon our inventory. For an accurate list of products in the current batch, please contact techserv@caymanchem.com.

### Contents of Bromodomain Inhibitor Set

Item No.	Item Name	Target(s)	Size
21033	ABBV-075	BRD2, BRD3, BRD4, BRDT	1 mg
20864	AZD 5153	BRD4	1 mg
19777	BAY-299	BRD1, TAF1	1 mg
22701	BAY-364	Negative control	1 mg
17448	BAZ2-ICR	BAZ2A, BAZ2B	1 mg
20311	BI-7273	BRD7, BRD9	1 mg
17897	BI-9564	BRD7, BRD9, CECR2	1 mg
14119	Bromosporine	Non-selective bromodomain inhibitor	1 mg
19956	CeMMEC1	TAF1	1 mg
20224	CeMMEC13	TAF1	1 mg
15479	CPI-203	BRD4	1 mg
23500	CPI-637	BRD9, CBP, p300	1 mg
14120	GSK2801	BAZ2A, BAZ2B	1 mg
18123	GSK5959	BRPF1	1 mg
20985	GSK6853	BRPF1	1 mg
11181	I-BET151	BRD2, BRD3, BRD4	1 mg
16872	I-BET726	BRD2, BRD3, BRD4	1 mg
17749	I-BRD9	BRD9	1 mg
14468	I-CBP112 (hydrochloride)	CBP, p300	1 mg
11232	(-)-JQ1	Negative control	1 mg
11187	(+)-JQ1	BRD4	1 mg
9002910	(+)-JQ1 (free acid)	BRD4	1 mg
22401	Lin28 1632	BRD4, CBP	1 mg
17661	LP99	BRD7, BRD9	1 mg
17662	NI-57	BRPF1B, BRPF2, BRPF3	1 mg
17124	OF-1	BRPF1B, BRPF2, BRPF3	1 mg

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/02/2023

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION

Page 2 of 2



Item No.	Item Name	Target(s)	Size
15947	OTX015	BRD2, BRD3, BRD4	1 mg
18811	PF-CBP1	CBP, p300	1 mg
11155	PFI-1	CBP, p300	1 mg
15267	PFI-3	PB1, SMARCA2, SMARCA4	1 mg
17663	PFI-4	BRPF1	1 mg
14469	SGC-CBP30	CBP, p300	1 mg
20030	TP-472	BRD7, BRD9	1 mg
17123	UMB-32	BRD4, TAF1, TAF1L	1 mg
19173	XMD8-92	BRD4	1 mg

## CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM