

# PRODUCT INFORMATION

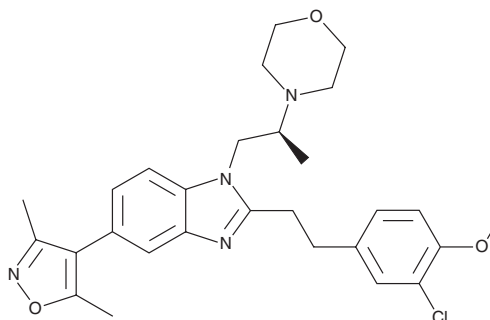


## SGC-CBP30

Item No. 14469

**CAS Registry No.:** 1613695-14-9  
**Formal Name:** 2-[2-(3-chloro-4-methoxyphenyl)ethyl]-5-(3,5-dimethyl-4-isoxazolyl)-1-[(2S)-2-(4-morpholinyl)propyl]-1H-benzimidazole

**MF:** C<sub>28</sub>H<sub>33</sub>ClN<sub>4</sub>O<sub>3</sub>  
**FW:** 509.0  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 221, 287 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

SGC-CBP30 is supplied as a crystalline solid. A stock solution may be made by dissolving the SGC-CBP30 in the solvent of choice, which should be purged with an inert gas. SGC-CBP30 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SGC-CBP30 in these solvents is approximately 5, 10, and 15 mg/ml, respectively.

SGC-CBP30 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SGC-CBP30 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SGC-CBP30 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

cAMP-responsive element-binding protein binding protein (CREBBP) and E1A-associated protein p300 (EP300) are transcriptional co-activators. Each contains a bromodomain, which recognizes acetylated lysine residues on target proteins. SGC-CBP30 is a potent inhibitor of CREBBP and EP300 bromodomains (IC<sub>50</sub>s = 21-69 and 38 nM, respectively).<sup>1</sup> Developed by the Structural Genomics Consortium (SGC), this compound displays over 40-fold selectivity for CREBBP over BRD4(1) and is functional in cells.

### Reference

1. Gallenkamp, D., Gelato, K.A., Haendler, B., *et al.* Bromodomains and their pharmacological inhibitors. *ChemMedChem* **9**(3), 438-464 (2014).

#### 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, 10/04/2022

#### 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