

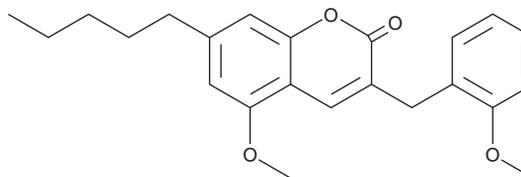
# PRODUCT INFORMATION



## PSB-SB1202

Item No. 15621

**CAS Registry No.:** 1399049-60-5  
**Formal Name:** 5-methoxy-3-[[2-methoxyphenyl)methyl]-7-pentyl-2H-1-benzopyran-2-one  
**MF:** C<sub>23</sub>H<sub>26</sub>O<sub>4</sub>  
**FW:** 366.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 251, 309 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

PSB-SB1202 is supplied as a crystalline solid. A stock solution may be made by dissolving the PSB-SB1202 in the solvent of choice. PSB-SB1202 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of PSB-SB1202 in ethanol and DMF is approximately 25 mg/ml and approximately 10 mg/ml in DMSO.

### Description

PSB-SB1202 is a benzylcoumarin compound that acts as an agonist at the major cannabinoid (CB) receptors, with EC<sub>50</sub> values of 56 and 14 nM at CB<sub>1</sub> and CB<sub>2</sub>, respectively.<sup>1</sup> It is highly selective for the CB<sub>1</sub> and CB<sub>2</sub> receptors (K<sub>i</sub>s = 32 and 49 nM, respectively), having minimal activity at GPR55 (IC<sub>50</sub> = 6.4 μM), a receptor that binds CBs and lysophosphatidylinositol.<sup>2</sup>

### References

1. Rempel, V., Volz, N., Hinz, S., *et al.* 7-Alkyl-3-benzylcoumarins: A versatile scaffold for the development of potent and selective cannabinoid receptor agonists and antagonists. *J. Med. Chem.* **55**(18), 7967-7977 (2012).
2. Rempel, V., Volz, N., Gläser, F., *et al.* Antagonists for the orphan G-protein-coupled receptor GPR55 based on a coumarin scaffold. *J. Med. Chem.* **56**(11), 4798-4810 (2013).

#### 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, 02/06/2024

#### 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