

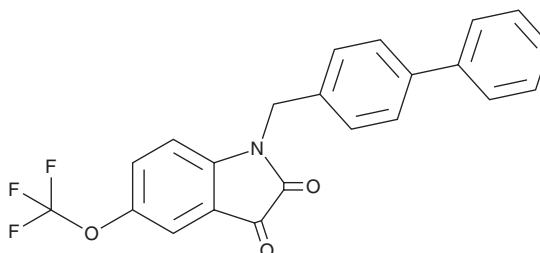
# PRODUCT INFORMATION



**VU0365114**

Item No. 33709

**CAS Registry No.:** 1208222-39-2  
**Formal Name:** 1-([1,1'-biphenyl]-4-ylmethyl)-5-(trifluoromethoxy)-1H-indole-2,3-dione  
**MF:** C<sub>22</sub>H<sub>14</sub>F<sub>3</sub>NO<sub>3</sub>  
**FW:** 397.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 248 nm  
**Supplied as:** A 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

VU0365114 is supplied as a solid. A stock solution may be made by dissolving the VU0365114 in the solvent of choice, which should be purged with an inert gas. VU0365114 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of VU0365114 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

## Description

VU0365114 is a positive allosteric modulator of M<sub>5</sub> muscarinic acetylcholine receptors (mAChRs).<sup>1</sup> It increases ACh-induced calcium mobilization in CHO cells expressing human M<sub>5</sub> mAChRs (EC<sub>50</sub> = 2.7 μM) selectively over CHO cells expressing human M<sub>1-4</sub> mAChRs (EC<sub>50</sub>s = >30 μM for all).

## Reference

1. Bridges, T.M., Kennedy, J.P., Hopkins, C.R., *et al.* Heterobiaryl and heterobiaryl ether derived M<sub>5</sub> positive allosteric modulators. *Bioorg. Med. Chem. Lett.* **20(19)**, 5617-5622 (2010).

### 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, 11/03/2022

## 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](http://WWW.CAYMANCHEM.COM)