

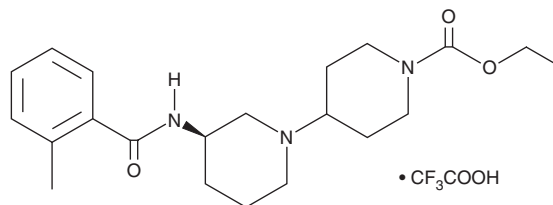
# PRODUCT INFORMATION



## VU0364572 (trifluoroacetate salt)

Item No. 20466

**CAS Registry No.:** 1240514-89-9  
**Formal Name:** (3R)-3-[(2-methylbenzoyl)amino]-[1,4'-bipiperidine]-1'-carboxylic acid, ethyl ester, 2,2,2-trifluoroacetate  
**MF:** C<sub>21</sub>H<sub>31</sub>N<sub>3</sub>O<sub>3</sub> • CF<sub>3</sub>COOH  
**FW:** 487.5  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Special Conditions:** Hygroscopic



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

### Laboratory Procedures

VU0364572 (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the VU0364572 (trifluoroacetate salt) in the solvent of choice. VU0364572 (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of VU0364572 (trifluoroacetate salt) in these solvents is approximately 25, 14, and 16 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of VU0364572 (trifluoroacetate salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of VU0364572 (trifluoroacetate salt) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

VU0364572 is a potent and selective agonist of the muscarinic acetylcholine receptor 1 (M<sub>1</sub>; EC<sub>50</sub> = 0.2 μM compared to >10 μM for M<sub>2</sub>-M<sub>5</sub>).<sup>1</sup> It induces concentration-dependent increases in Ca<sup>2+</sup> mobilization and ERK1/2 phosphorylation in CHO cells transfected with the human M<sub>1</sub> receptor (EC<sub>50</sub>s = 0.89 and 10 μM, respectively) with no effect on β-arrestin recruitment. VU0364572 also improves hippocampal-dependent spatial learning and localization abilities and acquisition of contextual fear learning but cannot reverse amphetamine-induced hyperlocomotion in rats. This suggests therapeutic potential of VU0364572 for the treatment of Alzheimer's disease but not psychosis.<sup>2</sup>

### References

1. Lebois, E.P., Digby, G.J., Sheffler, D.J., *et al.* Development of a highly selective, orally bioavailable and CNS penetrant M<sub>1</sub> agonist derived from the MLPCN probe ML071. *Bioorg. Med. Chem. Lett.* **21**(21), 6451-6455 (2011).
2. Digby, G.J., Noetzel, M.J., Bubser, M., *et al.* Novel allosteric agonists of M<sub>1</sub> muscarinic acetylcholine receptors induce brain region-specific responses that correspond with behavioral effects in animal models. *J. Neurosci.* **32**(25), 8532-8544 (2012).

#### 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, 06/05/2023

#### 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