

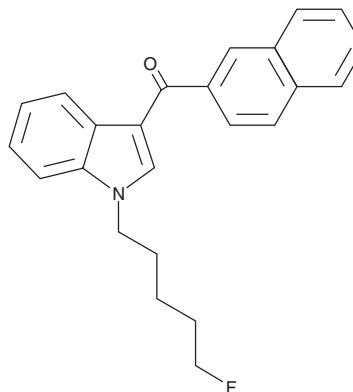
# PRODUCT INFORMATION



## AM2201 2'-naphthyl isomer

Item No. 10862

**CAS Registry No.:** 2365471-27-6  
**Formal Name:** (1-(5-fluoropentyl)-1H-indol-3-yl)  
(naphthalen-2-yl)methanone  
**MF:** C<sub>24</sub>H<sub>22</sub>FNO  
**FW:** 359.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 216, 252, 327 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Laboratory Procedures

AM 2201 2'-naphthyl isomer is supplied as a crystalline solid. A stock solution may be made by dissolving the AM2201 2'-naphthyl isomer in the solvent of choice. AM2201 2'-naphthyl isomer is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of AM2201 2'-naphthyl isomer in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

### Description

AM2201 is a potent synthetic cannabinoid (CB) with K<sub>i</sub> values of 1.0 and 2.6 nM for the CB<sub>1</sub> and CB<sub>2</sub> receptors, respectively.<sup>1</sup> AM2201 2'-naphthyl isomer differs structurally from AM2201 by having the naphthyl group attached at the 2' position. The biological and toxicological properties of this compound have not been characterized. This product is intended for forensic purposes.

### Reference

1. Makriyannis, A. and Deng, H. Cannabimimetic indole derivatives. WO 01/28557 A1 (2001), 1-25, PCT/US00/28832.

#### 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, 06/16/2022

#### 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