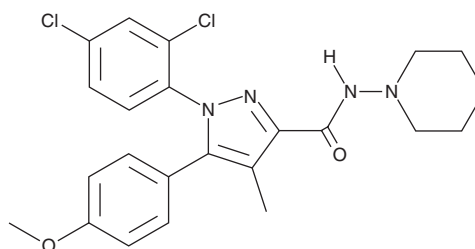


# PRODUCT INFORMATION

**NIDA-41020**

Item No. 21020

**CAS Registry No.:** 502486-89-7  
**Formal Name:** 1-(2,4-dichlorophenyl)-5-(4-methoxyphenyl)-4-methyl-N-1-piperidinyl-1H-pyrazole-3-carboxamide  
**MF:** C<sub>23</sub>H<sub>24</sub>Cl<sub>2</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 459.4  
**Purity:** ≥95%  
**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

NIDA-41020 is supplied as a crystalline solid. A stock solution may be made by dissolving the NIDA-41020 in the solvent of choice, which should be purged with an inert gas. NIDA-41020 is soluble in organic solvents such as ethanol and DMSO. The solubility of NIDA-41020 in these solvents is approximately 100 mM.

## Description

NIDA-41020 is a cannabinoid (CB) receptor 1 antagonist ( $K_i = 4.1$  nM).<sup>1</sup> It binds selectively to CB<sub>1</sub> over CB<sub>2</sub> ( $K_{bs} = 26$  and 831 nM, respectively).<sup>2</sup> NIDA-41020 was designed as a potential radioligand for use in positron emission tomography (PET).<sup>1</sup>

## References

1. Katoch-Rouse, R., Pavlova, O.A., Caulder, T., *et al.* Synthesis, structure-activity relationship, and evaluation of SR141716 analogues: Development of central cannabinoid receptor ligands with lower lipophilicity. *J. Med Chem.* **46**(4), 642-645 (2003).
2. Donohue, S.R., Halldin, C., and Pike, V.W. Synthesis and structure-activity relationships (SARs) of 1,5-diarylpyrazole cannabinoid type-1 (CB<sub>1</sub>) receptor ligands for potential use in molecular imaging. *Bioorg. Med. Chem.* **14**(11), 3712-3720 (2006).

### 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, 11/15/2022

## 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