

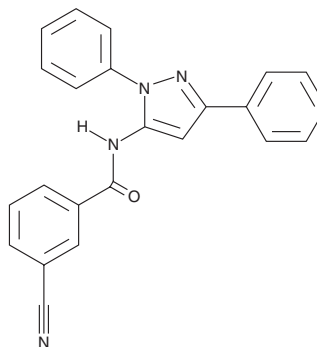
PRODUCT INFORMATION



CDPPB

Item No. 34132

CAS Registry No.: 781652-57-1
Formal Name: 3-cyano-N-(1,3-diphenyl-1H-pyrazol-5-yl)-benzamide
MF: C₂₃H₁₆N₄O
FW: 364.4
Purity: ≥98%
UV/Vis.: λ_{max}: 264 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

CDPPB is supplied as a solid. A stock solution may be made by dissolving the CDPPB in the solvent of choice, which should be purged with an inert gas. CDPPB is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of CDPPB in DMSO and DMF is approximately 25 and 20 mg/ml, respectively.

CDPPB is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CDPPB should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CDPPB has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CDPPB is a positive allosteric modulator (PAM) of metabotropic glutamate receptor 5 (mGluR5).¹ It potentiates glutamate-induced increases in intracellular calcium in CHO cells expressing mGluR5 (EC₅₀s = 10 and 20 nM for the human and rat receptor, respectively). CDPPB (30 mg/kg, s.c.) inhibits amphetamine-induced disruptions in prepulse inhibition (PPI) of the acoustic startle response in rats.

Reference

1. Lindsley, C.W., Wisnoski, D.D., Leister, W.H., *et al.* Discovery of positive allosteric modulators for the metabotropic glutamate receptor subtype 5 from a series of N-(1,3-diphenyl-1H-pyrazol-5-yl)benzamides that potentiate receptor function in vivo. *J. Med. Chem.* **47**(24), 5825-5828 (2004).

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, 10/25/2022

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