

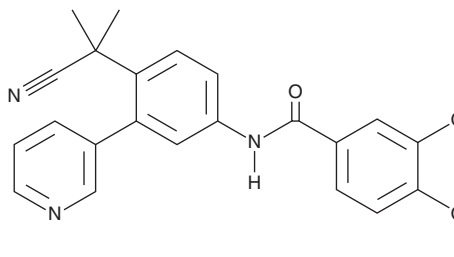
PRODUCT INFORMATION



ADX-68692

Item No. 46306

CAS Registry No.: 1067191-08-5
Formal Name: N-[4-(1-cyano-1-methylethyl)-3-(3-pyridinyl)phenyl]-3,4-dimethoxy-benzamide
MF: C₂₄H₂₃N₃O₃
FW: 401.5
Purity: ≥95%
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

ADX-68692 is supplied as a solid. A stock solution may be made by dissolving the ADX-68692 in the solvent of choice, which should be purged with an inert gas. ADX-68692 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

ADX-68692 is a negative allosteric modulator of the follicle-stimulating hormone receptor (FSHR) that reduces FSH-stimulated increases in progesterone and estradiol levels in primary rat granulosa cells (IC₅₀s = 1 and 0.82 μM, respectively).¹ It inhibits estrus and FSH-stimulated ovulation in female rats.

Reference

1. Dias, J.A., Campo, B., Weaver, B.A., *et al.* Inhibition of follicle-stimulating hormone-induced preovulatory follicles in rats treated with a nonsteroidal negative allosteric modulator of follicle-stimulating hormone receptor. *Biol. Reprod.* **90**(1), 19 (2014).

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, 07/31/2026

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