

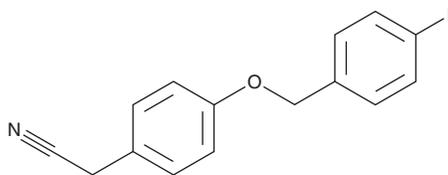
PRODUCT INFORMATION



MDK35833

Item No. 34332

CAS Registry No.: 1016535-83-3
Formal Name: 4-[(4-fluorophenyl)methoxy]-benzeneacetonitrile
MF: C₁₅H₁₂FNO
FW: 241.3
Purity: ≥98%
UV/Vis.: λ_{max}: 227 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

MDK35833 is supplied as a solid. A stock solution may be made by dissolving the MDK35833 in the solvent of choice, which should be purged with an inert gas. MDK35833 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MDK35833 in ethanol is approximately 15 mg/ml and approximately 30 mg/ml in DMSO and DMF.

MDK35833 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MDK35833 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MDK35833 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MDK35833 is an inducer of octamer-binding transcription factor 3/4 (Oct3/4) expression.¹ It induces Oct3/4 expression in reporter assays using HEK or NCCIT cells when used at a concentration of 10 μM. MDK35833 (20 μM) induces expression of the genes encoding the pluripotency-associated proteins NANOG, REX1, and KLF4 in primary human fibroblasts.

Reference

1. Cheng, X., Dimou, E., Alborzinia, H., *et al.* Identification of 2-[4-[(4-Methoxyphenyl)methoxy]-phenyl] acetonitrile and derivatives as potent Oct3/4 inducers. *J. Med. Chem.* **58**(12), 4976-4983 (2015).

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, 11/22/2022

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