

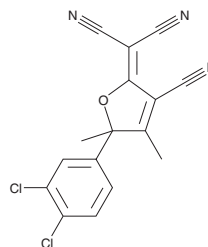
PRODUCT INFORMATION



YE 120

Item No. 29720

CAS Registry No.: 383124-82-1
Formal Name: 2-[3-cyano-5-(3,4-dichlorophenyl)-4,5-dimethyl-2(5H)-furanylidene]-propanedinitrile
MF: C₁₆H₉Cl₂N₃O
FW: 330.2
Purity: ≥98%
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

YE 120 is supplied as a solid. A stock solution may be made by dissolving the YE 120 in the solvent of choice, which should be purged with an inert gas. YE 120 is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

YE 120 is an agonist of G protein-coupled receptor 35 (GPR35) with an EC₅₀ value of 32 nM in a dynamic mass redistribution (DMR) assay using HT-29 cells.¹ It induces β-arrestin translocation in Tango GPR35-*bla* U2OS cells (EC₅₀ = 10.2 μM). YE 120 (10 μM) promotes wound closure in YAMC colonic epithelial and IEC-6 intestinal epithelial cell monolayers, an effect that can be blocked by the GPR35 antagonist CID-2745687 (Item No. 12046).²

References

1. Deng, H., Hu, H., He, M., *et al.* Discovery of 2-(4-methylfuran-2(5H)-ylidene)malononitrile and thieno[3,2-b]thiophene-2-carboxylic acid derivatives as G protein-coupled receptor 35 (GPR35) agonists. *J. Med. Chem.* **54**(20), 7385-7396 (2011).
2. Tsukahara, T., Hamouda, N., Utsumi, D., *et al.* G protein-coupled receptor 35 contributes to mucosal repair in mice via migration of colonic epithelial cells. *Pharmacol. Res.* **123**, 27-39 (2017).

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, 12/08/2022

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