

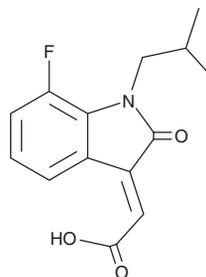
PRODUCT INFORMATION



ASP7663

Item No. 31122

CAS Registry No.: 1190217-35-6
Formal Name: (2E)-2-[7-fluoro-1,2-dihydro-1-(2-methylpropyl)-2-oxo-3H-indol-3-ylidene]-acetic acid
MF: C₁₄H₁₄FNO₃
FW: 263.3
Purity: ≥95%
UV/Vis.: λ_{max}: 224, 250, 255, 315 nm
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

ASP7663 is supplied as a crystalline solid. A stock solution may be made by dissolving the ASP7663 in the solvent of choice, which should be purged with an inert gas. ASP7663 is soluble in organic solvents such as ethanol and DMSO. The solubility of ASP7663 in these solvents is approximately 25 mg/ml.

Description

ASP7663 is a transient receptor potential ankyrin 1 (TRPA1) agonist.¹ It induces calcium influx in HEK293 cells expressing human, rat, or mouse TRPA1 (EC₅₀s = 0.51, 0.54, and 0.5 μM, respectively) and stimulates serotonin (5-HT) release from QGP-1 enterochromaffin cells that endogenously express TRPA1 (EC₅₀ = 72.5 μM). It reduces colonic transit time in mouse models of constipation induced by clonidine (Item No. 15949) or loperamide when administered orally at a dose of 1 mg/kg, an effect that can be blocked by vagotomy or the TRPA1 antagonist HC-030031 (Item No. 11923). ASP7663 (0.3 mg/kg, p.o.) reduces the number of abdominal contractions induced by colorectal distension in mice.

Reference

1. Kojima, R., Nozawa, K., Doihara, H., *et al.* Effects of novel TRPA1 receptor agonist ASP7663 in models of drug-induced constipation and visceral pain. *Eur. J. Pharmacol.* **723**, 288-293 (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, 12/12/2022

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