

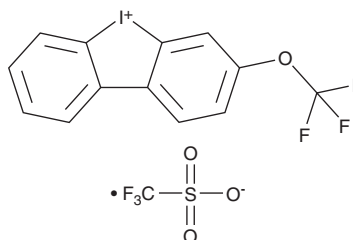
PRODUCT INFORMATION



GPR3 Agonist 2

Item No. 42616

CAS Registry No.: 1983120-51-9
Formal Name: 3-(trifluoromethoxy)-dibenziodolium,
1,1,1-trifluoromethanesulfonate
Synonym: G Protein-Coupled Receptor 3 Agonist 2
MF: $C_{13}H_7F_3IO \cdot CF_3O_3S$
FW: 512.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GPR3 agonist 2 is supplied as a solid. A stock solution may be made by dissolving the GPR3 agonist 2 in the solvent of choice, which should be purged with an inert gas. GPR3 agonist 2 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

GPR3 agonist 2 is an agonist of G protein-coupled receptor 3 (GPR3).¹ It induces cAMP accumulation in HEK293 cells expressing human GPR3 ($EC_{50} = 0.26 \mu M$). It is selective for GPR3 over GPR6 and GPR12 at $10 \mu M$.

Reference

1. Gay, E.A., Harris, D.L., Wilson, J.W., *et al.* The development of diphenyleneiodonium analogs as GPR3 agonists. *Bioorg. Med. Chem. Lett.* **94**, 129427 (2023).

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, 02/27/2025

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