

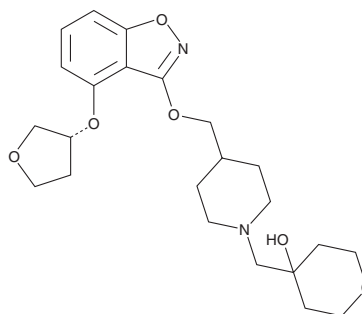
PRODUCT INFORMATION



PF-04995274

Item No. 28776

CAS Registry No.: 1331782-27-4
Formal Name: tetrahydro-4-[[[4-[[[(3R)-tetrahydro-3-furanyl]oxy]-1,2-benzisoxazol-3-yl]oxy]methyl]-1-piperidinyl]methyl]-2H-pyran-4-ol
MF: C₂₃H₃₂N₂O₆
FW: 432.5
Purity: ≥98%
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

PF-04995274 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-04995274 in the solvent of choice, which should be purged with an inert gas. PF-04995274 is soluble in organic solvents such as chloroform and dimethyl formamide (DMF). The solubility of PF-04995274 in these solvents is approximately 30 and 20 mg/ml, respectively.

PF-04995274 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PF-04995274 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PF-04995274 has a solubility of approximately 0.25 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PF-04995274 is a partial agonist of the serotonin (5-HT) receptor subtypes 5-HT_{4A}, 5-HT_{4B}, 5-HT_{4D}, and 5-HT_{4E} (K_is = 0.36, 0.46, 0.15, and 0.32 nM, respectively, in a radioligand binding assay).¹ It increases cAMP levels in HEK293 cells expressing 5-HT_{4A}, 5-HT_{4B}, 5-HT_{4D}, and 5-HT_{4E} receptors (EC₅₀s = 0.47, 0.36, 0.37, and 0.26 nM, respectively, for the human recombinant receptors). PF-04995274 decreases scopolamine-induced increases in the distance traveled in the Morris water maze in rats when administered at a dose of 0.032 mg/kg.

Reference

1. Brodney, M. A., Johnson, D. E., Sawant-Basak, A. *et. al.* Identification of multiple 5-HT₄ partial agonist clinical candidates for the treatment of alzheimer's disease. *J. Med. Chem.* **55**, 9240-9254 (2012).

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, 12/06/2022

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