

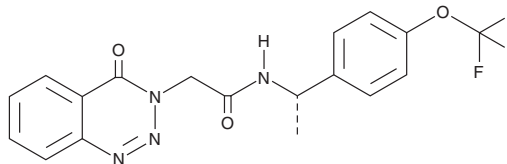
PRODUCT INFORMATION



TAK-041

Item No. 36108

CAS Registry No.: 1929519-13-0
Formal Name: 4-oxo-N-[(1S)-1-[4-(trifluoromethoxy)phenyl]ethyl]-1,2,3-benzotriazine-3(4H)-acetamide
Synonym: NBI 1065864
MF: C₁₈H₁₅F₃N₄O₃
FW: 392.3
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

TAK-041 is supplied as a crystalline solid. A stock solution may be made by dissolving the TAK-041 in the solvent of choice, which should be purged with an inert gas. TAK-041 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of TAK-041 in these solvents is approximately 1 and 3 mg/ml, respectively.

Description

TAK-041 is an agonist of G protein-coupled receptor 139 (GPR139).¹ It induces calcium mobilization in cells expressing human GPR139 (EC₅₀ = 22 nM). *In vivo*, TAK-041 (0.3 and 3 mg/kg) increases social interaction in mice.

Reference

1. Reichard, H.A., Schiffer, H.H., Monenschein, H., *et al.* Discovery of TAK-041: A potent and selective GPR139 agonist explored for the treatment of negative symptoms associated with schizophrenia. *J. Med. Chem.* **64**(15), 11527-11542 (2021).

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, 11/08/2022

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