

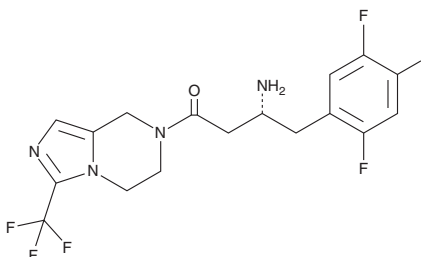
PRODUCT INFORMATION



CGT-8012

Item No. 45055

CAS Registry No.: 1152810-23-5
Formal Name: 3R-amino-1-[5,6-dihydro-3-(trifluoromethyl)imidazo[1,5-a]pyrazin-7(8H)-yl]-4-(2,4,5-trifluorophenyl)-1-butanone
Synonym: Sheng Gelieting
MF: C₁₇H₁₆F₆N₄O
FW: 406.3
Purity: ≥95%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

CGT-8012 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and acetonitrile purged with an inert gas can be used. CGT-8012 is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in acetonitrile.

Description

CGT-8012 is a dipeptidyl peptidase 4 (DPP-4) inhibitor (IC₅₀ = 0.14 μM).¹

Reference

1. Tang, P.C., Yang, F., Wang, Y., et al. Piperazine derivatives, preparation process and pharmaceutical use thereof. *Shanghai Hengrui Pharmaceutical Co. Ltd. WO 2009/065298 A1* (2007).

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, 03/10/2026

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