

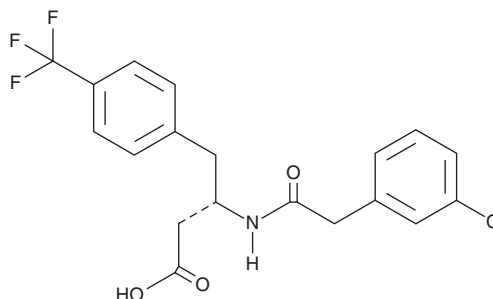
PRODUCT INFORMATION



CATPB

Item No. 45837

CAS Registry No.: 1322598-09-3
Formal Name: β S-[[2-(3-chlorophenyl)acetyl]amino]-4(trifluoromethyl)-benzenebutanoic acid
MF: $C_{19}H_{17}ClF_3NO_3$
FW: 399.8
Purity: $\geq 98\%$
Supplied as: A 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

CATPB is supplied as a solid. A stock solution may be made by dissolving the CATPB in the solvent of choice, which should be purged with an inert gas. CATPB is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

CATPB is a free fatty acid receptor 2 (FFAR2/GPR43) antagonist.¹ It inhibits agonist-induced activation of FFAR2/GPR43 in cells expressing FFAR2/GPR43 ($IC_{50} = 224$ nM). CATPB (500 nM) reduces agonist-induced superoxide production in neutrophils primed with TNF- α or latrunculin A (Item No. 10010630).²

References

1. Højgaard Hansen, A., Christensen, H.B., Pandey, S.K., *et al.* Structure-activity relationship explorations and discovery of a potent antagonist for the free fatty acid receptor 2. *ChemMedChem* **16**(21), 3326-3341 (2021).
2. Björkman, L., Mårtensson, J., Winther, M., *et al.* The neutrophil response induced by an agonist for free fatty acid receptor 2 (GPR43) is primed by tumor necrosis factor alpha and by receptor uncoupling from the cytoskeleton but attenuated by tissue recruitment. *Mol. Cell Biol.* **36**(20), 2583-2595 (2016).

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, 06/17/2026

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