

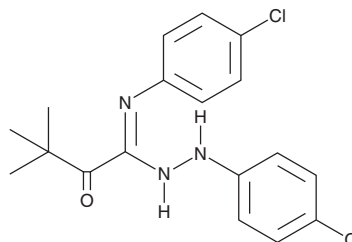
PRODUCT INFORMATION



TY 52156

Item No. 19119

CAS Registry No.: 934369-14-9
Formal Name: N-(4-chlorophenyl)-3,3-dimethyl-2-oxobutanimidic acid, 2-(4-chlorophenyl)hydrazide
MF: C₁₈H₁₉Cl₂N₃O
FW: 364.3
Purity: ≥98%
UV/Vis.: λ_{max}: 248, 301, 359 nm
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

TY 52156 is supplied as a crystalline solid. A stock solution may be made by dissolving the TY 52156 in the solvent of choice. TY 52156 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of TY 52156 in these solvents is approximately 15, 25, and 30 mg/ml, respectively.

TY 52156 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

TY 52156 is a selective sphingosine-1-phosphate receptor 3 (S1P₃) antagonist (K_i = 110 nM).¹ It preferentially inhibits the S1P-induced increase in intracellular calcium in Chinese hamster ovary cells expressing S1P₃ over cells expressing S1P₁, S1P₂, or S1P₄.¹ TY 52156 has been shown to suppress FTY720-induced S1P₃ receptor-mediated coronary flow in isolated perfused rat hearts.¹ TY 52156 can also inhibit S1P-induced breast cancer stem cell expansion *in vitro*.²

References

1. Murakami, A., Takasugi, H., Ohnuma, S., *et al.* Sphingosine 1-phosphate (S1P) regulates vascular contraction via S1P₃ receptor: Investigation based on a new S1P₃ receptor antagonist. *Mol. Pharmacol.* **77**(4), 704-713 (2016).
2. Hirata, N., Yamada, S., Shoda, T., *et al.* Sphingosine-1-phosphate promotes expansion of cancer stem cells via S1P₃ by a ligand-independent Notch activation. *Nat. Commun.* **5**:4806, (2014).

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, 10/31/2022

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