PRODUCT INFORMATION



GR148672X

Item No. 28756

CAS Registry No.:	263890-70-6	\sim
Formal Name:	4,4,4-trifluoro-1-(2-thienyl)-1,2,3-butanetrione,	
	2-[(3-methylphenyl)hydrazone]	
MF:	$C_{15}H_{11}F_{3}N_{2}O_{2}S$	F N N
FW:	340.3	F L
Purity:	≥98%	F H
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

GR148672X is supplied as a solid. A stock solution may be made by dissolving the GR148672X in the solvent of choice, which should be purged with an inert gas. GR148672X is soluble in DMSO.

Description

GR148672X is an inhibitor of triacylglycerol hydrolase/carboxylesterase 1 (TGH/CES1; IC₅₀ = 4 nM for the human hepatic enzyme).^{1,2} It is selective for TGH/CES1 over lipoprotein lipase (LPL) at 5 $\mu M.^1$ GR148672X (25 mg/kg twice per day) decreases plasma levels of triglycerides, total, VLDL, and LDL cholesterol, and apolipoprotein B-100 (ApoB-100) in hamsters.

References

- 1. Borg-Capra, C.S., Lehner, R.J., and Vance, D.E. Method of screening for triacyglycerol hydrolase inhibitors. Glaxo Group Limited and Governors of the University of Alberta. WO 01/16358 A2 (2001).
- 2. Wei, E., Ben Ali, Y., Lyon, J., et al. Loss of TGH/Ces3 in mice decreases blood lipids, improves glucose tolerance, and increases energy expenditure. Cell Metab. 11(3), 183-193 (2010).

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.

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, 11/03/2022

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