

PRODUCT INFORMATION



GW 501516

Item No. 36516

CAS Registry No.: 317318-70-0

Formal Name: 2-[2-methyl-4-[[[4-methyl-2-[4-(trifluoromethyl)phenyl]-5-thiazolyl]methyl]thio]phenoxy]-acetic acid

Synonym: GW 1516

MF: C₂₁H₁₈F₃NO₃S₂

FW: 453.5

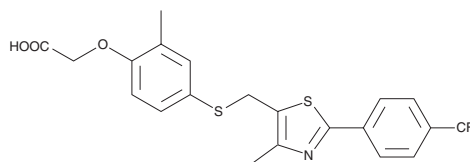
Purity: ≥98%

UV/Vis.: λ_{max}: 318 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥5 years



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

Description

GW 501516 (Item No. 36516) is an analytical reference standard categorized as a peroxisome proliferator-activated receptor δ (PPAR δ) agonist.¹ Formulations containing GW 501516 have been used to enhance physical performance in athletes. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 10004272).

Reference

1. Thevis, M., Möller, I., Thomas, A., *et al.* Characterization of two major urinary metabolites of the PPAR δ -agonist GW1516 and implementation of the drug in routine doping controls. *Anal. Bioanal. Chem.* **396**(7), 2479-2491 (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.

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, 07/07/2023

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