

PRODUCT INFORMATION



GW 6471

Item No. 11697

CAS Registry No.: 880635-03-0

Formal Name: N-[(2S)-2-[[[(1Z)-1-methyl-3-oxo-3-[4-(trifluoromethyl)phenyl]-1-propen-1-yl]amino]-3-[4-[2-(5-methyl-2-phenyl-4-oxazolyl)ethoxy]phenyl]propyl]-propanamide

MF: $C_{35}H_{36}F_3N_3O_4$

FW: 619.7

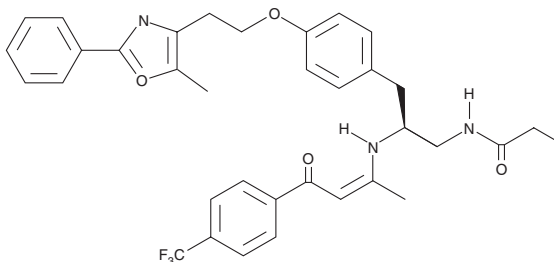
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 225, 278, 347, 352 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

GW 6471 is supplied as a crystalline solid. A stock solution may be made by dissolving the GW 6471 in the solvent of choice, which should be purged with an inert gas. GW 6471 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GW 6471 in these solvents is approximately 10, 11, and 20 mg/ml, respectively.

GW 6471 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GW 6471 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. GW 6471 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The peroxisome proliferator-activated receptor α (PPAR α) is a nuclear receptor that regulates the expression of genes involved in fatty acid metabolism, lipoprotein synthesis and metabolism, and inflammation.¹⁻³ GW 6471 is an antagonist of PPAR α ($IC_{50} = 240$ nM).⁴ It drives the displacement of coactivators from PPAR α and promotes the recruitment of co-repressor proteins like nuclear co-repressor.⁴ GW 6471 is used in cell-free, cell lysate, and whole cell systems to examine the impact of PPAR α antagonism.⁵⁻⁷

References

1. Kersten, S., Desvergne, B., and Wahli, W. *Nature* **405**, 421-424 (2000).
2. Tenenbaum, A., Motro, M., and Fisman, E.Z. *Cardiovasc. Diabetol.* **4**(14), 1-5 (2005).
3. Ahmed, W., Ziouzenkova, O., Brown, J., et al. *J. Intern. Med.* **262**, 184-198 (2007).
4. Xu, H.E., Stanley, T.B., Montana, V.G., et al. *Nature* **415**(6873), 813-817 (2002).
5. Wang, W., Lin, Q., Lin, R., et al. *Exp. Cell Res.* **319**(10), 1523-1533 (2013).
6. Downer, E.J., Clifford, E., Amu, S., et al. *J. Biol. Chem.* **287**(30), 25440-25453 (2012).
7. Müller, M.Q., de Koning, L.I., Schmidt, A., et al. *J. Med. Chem.* **52**(9), 2875-2879 (2009).

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, 09/29/2022

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