

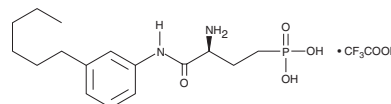
PRODUCT INFORMATION



W140 (trifluoroacetate salt)

Item No. 10009110

CAS Registry No.: 909725-64-0
Formal Name: [3S-amino-4-[(3-hexylphenyl)amino]-4-oxobutyl]-
phosphonic acid, mono(trifluoroacetate)
MF: C₁₆H₂₇N₂O₄P • CF₃COOH
FW: 456.4
Purity: ≥95%
UV/Vis.: λ_{max}: 207, 246 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

W140 (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the W140 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. W140 (trifluoroacetate salt) is soluble in the organic solvent ethanol at a concentration of approximately 0.15 mg/ml.

Description

W140 is the inactive enantiomer of the sphingosine-1-phosphate receptor 1 (S1P₁) antagonist W146 (Item No. 10009109).¹ It only weakly binds to S1P₁ (K_i = 4.6 μM for the human receptor). W140 has been used as a negative control *in vitro* and *in vivo* in the study of S1P₁ signaling.^{1,2}

References

1. Sanna, M.G., Wang, S.K., Gonzalez-Cabrera, P.J., *et al.* Enhancement of capillary leakage and restoration of lymphocyte egress by a chiral S1P₁ antagonist *in vivo*. *Nat. Chem. Biol.* **2(8)**, 434-441 (2006).
2. Li, C., Chi, X.X., Xie, W., *et al.* Sphingosine 1-phosphate receptor 2 antagonist JTE-013 increases the excitability of sensory neurons independently of the receptor. *J. Neurophysiol.* **108(5)**, 1473-1483 (2012).

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, 01/25/2024

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