

AK-7

Epigenetic Enzyme Inhibitor

Catalog # S36-904

Lot # P3320-3

CAS # 420831-40-9

Product Description

Molecular Formula: C₁₉H₂₁BrN₂O₃S

Appearance: White powder

Melting Point: 148-151°C

Molecular Weight: 437.4

Purity: 98% (TLC); NMR (Conforms)

Solubilization: May be dissolved in DMSO (50 mg/ml); or ethanol (8 mg/ml, warm).

Alias

N-(3-Bromophenyl)-3-[(hexahydro-1H-azepin-1-yl)sulfonyl]benzamide

Specific Activity

SIRT2 Inhibitor

Storage and Stability

Store desiccated as supplied at ambient temperature for up to 3 years. Store solutions at -20°C for up to 3 months.

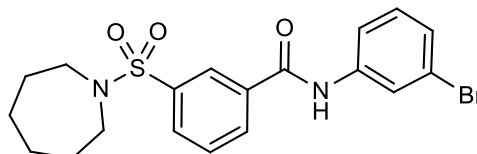
Scientific Background

A selective Sirt2 inhibitor (IC₅₀=15.5 μM) which does not inhibit Sirt1 or Sirt3¹. It down-regulates cholesterol biosynthesis genes in primary striatal neurons¹. Displays neuroprotective effects in Huntington's disease mouse models including improved motor function, extended survival and reduced brain atrophy all of which were associated with reduced aggregated mutant huntingtin².

References

1. DM Taylor et al. ACS Chem. Biol. 2011 6:540
2. V Chopra et al. Cell Rep. 2012 2:1492

Molecular Structure



AK-7

Epigenetic Enzyme Inhibitor

Catalog #	S36-904
CAS #	420831-40-9
Lot #	P3320-3
Purity	98% (TLC)
Format	White powder
Solubilization	May be dissolved in DMSO (50 mg/ml); or ethanol (8 mg/ml, warm)
Storage & Shipping	Store desiccated as supplied at ambient temperature for up to 3 years. Store solutions at -20°C for up to 3 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.