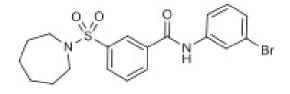


Catalog No: 14054, 14055 Format: 25 mg, 5 mg

Chemical Properties:

$$\begin{split} \mathsf{MW} &= 437.4\\ \mathsf{C}_{19}\mathsf{H}_{21}\mathsf{BrN}_2\mathsf{O}_3\mathsf{S}\\ \mathsf{CAS} & 420831\text{-}40\text{-}9\\ \mathsf{Physical Properties: White powder, MP = 148\text{-}151^\circ\mathsf{C}\\ \mathbf{Names:} \ \mathsf{AK}\text{-}7, \ 3\text{-}(1\text{-}Azepanylsulfonyl)\text{-}N\text{-}(3\text{-}bromophenyl)benzamide} \end{split}$$



Pharmacology: A selective Sirt2 inhibitor ($IC_{50} = 15.5$ M) which does not bind Sirt1 or Sirt3 (ref 1). It down-regulates cholesterol biosynthesis genes in primary striatal neurons (ref 1). Displays neuroprotective effects in Huntington's disease mouse models including improved motor function, extended survival and reduce brain atrophy all of which were associated with reduced aggregated mutant huntingtin (ref 2).

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

Fluorescent Properties: N/A

Quality Control: >98% (TLC); NMR (Conforms)

References:

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

Storage and Guarantee: Store desiccated as supplied at room temperature for up to 3 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.