

TC-E-5003

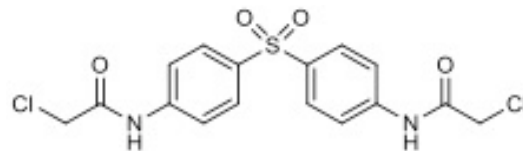
**Catalog No:** 14099**Format:** 50 mg**Chemical Properties:**

MW = 401.3

C<sub>16</sub>H<sub>14</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>4</sub>S

CAS 17328-16-4

Physical Properties: White powder

**Names:** N,N'-(Sulfonyldi-4,1-phenylene)bis(2-chloroacetamide)

**Pharmacology:** Selective protein arginine methyltransferase 1 (PRMT1) inhibitor (IC<sub>50</sub> = 1.5 μM). Displays no activity against CARM1 and Set7/9 methyltransferases. Inhibits growth of MCF7 breast cancer cells and LNCaP prostate cancer cells. Blocks androgen-induced gene expression in LNCaP cells (ref1).

**Solubilization:** May be dissolved in DMSO (50 mg/ml); DMF (50 mg/ml); Ethanol (nil)

**Fluorescent Properties:** N/A

**Quality Control:**

&gt;98% (TLC); NMR (Conforms)

## References:

1. EM Bissinger *et al. Bioorg. Med. Chem. Lett.* 2011, 19:3717

**Storage and Guarantee:** Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.