# **RS-1**

#### RAD51 activator / CRISPR enhancer

RS-1 is a cell-permeable sulfonamido-benzamide-based allosteric regulator. It enhances the DNA binding and recombinase activities of the human homologous recombination protein RAD51 by stabilizing its active conformation without affecting its ATP hydrolysis activity.

Key features and applications include:

- DNA Repair Studies: RS-1 promotes the formation of active RAD51 filaments on single-stranded DNA (ssDNA), enhancing homologous recombination efficiency.
- Genome Editing: It significantly increases the efficiency of CRISPR-Cas9 and TALEN-mediated genome editing by enhancing homologydirected repair (HDR).
- Cancer Research: RS-1 has been shown to increase the resistance of primary human neonatal dermal fibroblasts to cisplatin-induced cell death.

Relevant disease states include:

- **Cancer:** RS-1 has shown potential in promoting antitumor responses by enhancing DNA repair mechanisms, making it relevant for research into various cancers.
- **Genetic Disorders:** Its role in enhancing HDR makes it useful in studying genetic disorders where precise gene correction is needed.

## **Ordering Information**

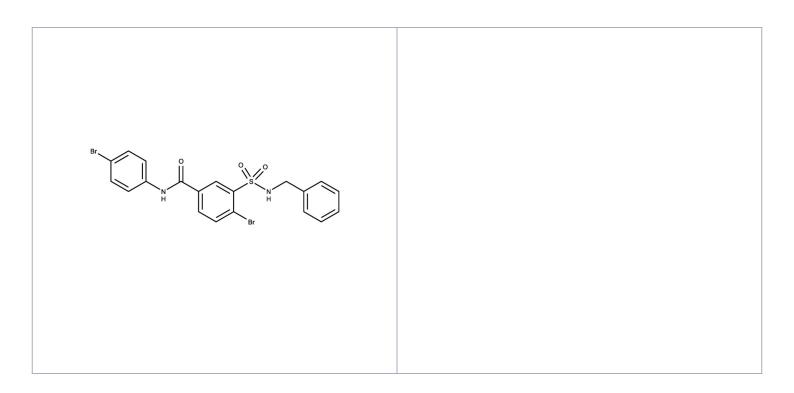
**Order Online** »

**ENZ-CHM337-0050** 50mg

Manuals, SDS & CofA

View Online »

· GMP format available



### **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended.

Solutions in DMSO may be stored at -20°C for up to 3 months.

**Handling** Keep container tightly closed in a dry and well-ventilated place.

Short Term Storage Ambient

Long Term Storage Ambient

**Shipping** Ambient Temperature

#### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name 4-Bromo-N-(4-bromophenyl)-3-

[[(phenylmethyl)amino]sulfonyl]benzamide

Appearance White solid.

CAS 312756-74-4

Couple Target Rad

Couple Type Activator

Formula  $C_{20}H_{16}Br_2N_2O_3S$ 

**Identity** Determined by NMR.

MW 524.23

**MeltingPoint** 200.8-202.4°C (393.4-396.3°F)

Purity ≥99% (HPLC)

**Solubility** Soluble in DMSO (30 mg/mL).



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989

eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com