

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 homology-directed 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

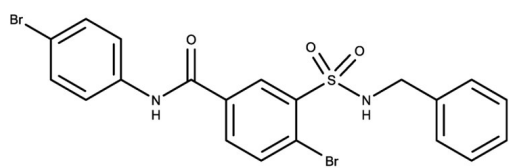
[Order Online »](#)

| | |
|-----------------|------|
| 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). |



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)