
DR2313

(Catalog No. M75002)

Alternative Names

2-Methyl-3,5,7,8-tetrahydrothiopyrano[4,3-d]pyrimidine-4-one

Description

PARP Inhibitor XI, DR2313 is a water soluble and brain-permeable pyrimidinone compound that functions as a Poly(ADP-ribose) polymerase inhibitor. Poly(ADP-ribose) polymerases, known as PARP, function by triggering DNA repair mechanisms and activating apoptosis in damaged cells. Studies show that DR2313 is a NAD⁺ utilizing enzyme which competitively inhibits Poly(ADP-ribose) polymerases but does not display selectivity among PARP-1 and PARP-2. In addition, in vitro studies of cerebral ischemia exhibiting an overactivation of PARP can be reduced by introducing DR2313

CAS Number

284028-90-6

Chemical Formula

C₈H₁₀N₂OS

Molecular Weight

182.2

Size

10 mg

Purity

>98%

Solubility

H₂O

Storage Temp

-20°C

Ordering Information:

| Products | Size | Cat. No. |
|----------|-------|----------|
| DR2313 | 10 mg | M75002-1 |

This product is for research purposes only. Not intended for use in diagnostic procedures.