

Geldanamycin, FITC-labeled

Catalog #: 50312

Lot #: -

Size: 30 µl

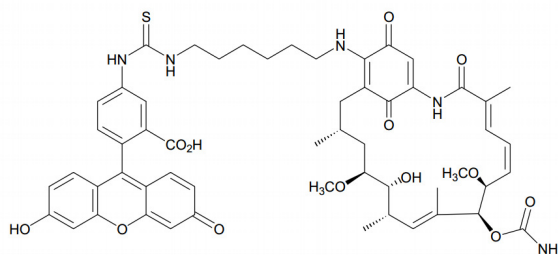
Structure:

CAS Registry #: 30562-34-6
(Geldanamycin)

Purity: ≥ 98%

Chemical Formula: C₅₅H₆₃N₅O₁₃S

Molecular Weight: 1034.2

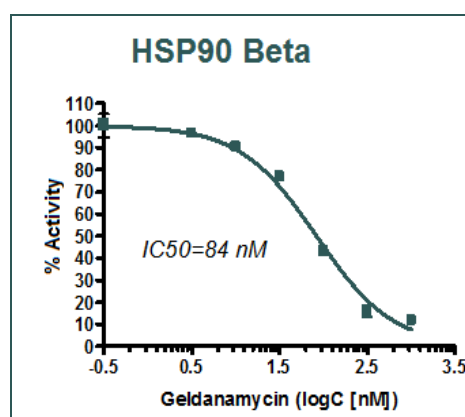
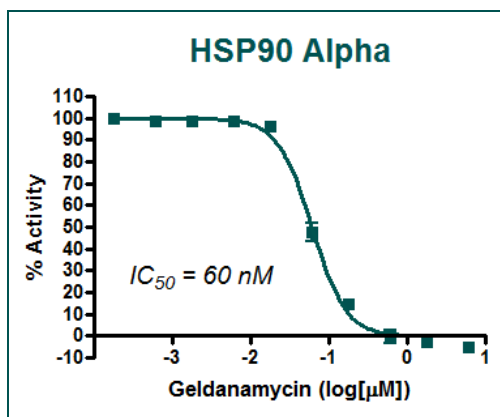


Description: A novel geldanamycin fluorescent probe optimized for use with BPS's HSP90 Assay Kits. May also be used for detection of cell surface HSP90.

Appearance: liquid

Supplied as: DMSO solution (2.5 µM)

Biological Activity: FITC-labeled geldanamycin has been shown to inhibit HSP90α protein with an IC₅₀ of 60 nM and HSP90β with an IC₅₀ of 84 nM as shown. Measured using the HSP90 Assay Kits.



Storage/Stability: >6 months at -80°C. Avoid freeze/thaw cycles. Keep DMSO stock solutions out of contact with air. Protect from light.

References:

1. Kim, J., *et al.*, *Biomol. Screening* 2004; **9(5)**: 375-381.
2. Howes, R., *et al.*, *Anal. Biochem.* 2006; **350**:202-213.