

17-NHS-ALA-GA

GA analogue for coupling of NH₂-containing molecules

Catalog # ant-nhg1-1

For research use only

Version # 16D07-MM

PRODUCT INFORMATION

Contents:

1 mg of 17-NHS-ALA-GA is supplied as a purple powder.

Storage and stability:

- 17-NHS-ALA-GA is shipped at room temperature. Store powder at -20 °C. Protect from light. Product as powder is stable for 6 months at -20 °C when properly stored.

- Once solubilized 17-NHS-ALA-GA should be aliquoted and stored at -20 °C. Protect from light. Resuspended product is stable 3 months at -20 °C when properly stored.

QUALITY CONTROL

- Purity: ≥95% (UHPLC).

DESCRIPTION

17-NHS-ALA-GA is an NHS (N-hydroxysuccinimide) activated geldanamycin analogue designed for easy coupling of NH₂-containing molecules. NHS coupling forms a chemically stable amide bond with ligands containing primary amino groups such as small proteins and peptides. Thus, 17-NHS-ALA-GA can be used for conjugation of geldanamycin to a monoclonal antibody similarly to 17-GMB-APA-GA^{1,2}. It can also be used to perform affinity chromatography to purify GA binding proteins such as Hsp90.

1. . Mandler R. et al., 2002. Modifications in synthesis strategy improve the yield and efficacy of geldanamycin-herceptin immunoconjugates. *Bioconj Chem.* 13(4):786-91. 2. Mandler R et al. 2000. Immunoconjugates of Geldanamycin and Anti-HER2 Monoclonal Antibodies.

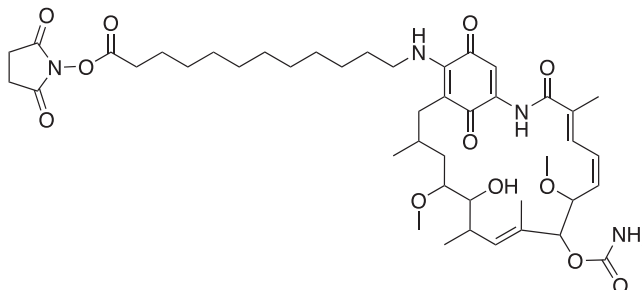
CHEMICAL PROPERTIES

Formula: C₄₄H₈₄N₄O₁₂

Molecular weight: 841

Solubility: DMSO (10 mg/ml)

Structure:



TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Hong Kong: +852 3-622-34-80
E-mail: info@invivogen.com