

## Product Information

## Description

CL2A-SN38 is a SN38 derivative with a Mal-PEG attached peptide linker, it can easily react with antibody to form an antibody-drug conjugate (ADC). SN-38 is a potent DNA Topoisomerase I inhibitor. CL2A-SN-38 provides significant and specific antitumor effects against a range of human solid tumor types. CL2A linker is cleavable through pH sensitivity, giving rise to bystander effect, and binds the antibody at a cysteine residue via a disulfide bond.

## Molecular Formula:

 $C_{73}H_{97}N_{11}O_{22}$ 

## Supplied As:

Solid

## Stability:

At least 6 months at -20°C. Avoid freeze/thaw cycles.

## Storage:

-20°C

## MW:

1480.6 Da

## Cas No.

1279680-68-0

## Chemical Structure

