

## Product Information

|                           |  |
|---------------------------|--|
| <b>Description</b>        | Auristatin PE is a novel synthetic Dolastatin 10 derivative. Auristatin PE inhibits tubulin polymerization, resulting in cell cycle arrest and induction of apoptosis. |
| <b>Molecular Formula:</b> | C <sub>39</sub> H <sub>67</sub> N <sub>5</sub> O <sub>6</sub>  |
| <b>Supplied As:</b>       | Solid  |
| <b>Stability:</b>         | At least 6 months at -20°C. Avoid freeze/thaw cycles.  |
| <b>Storage:</b>           | -20°C  |
| <b>MW:</b>                | 702.0 Da   |
| <b>Cas No.</b>            | 149606-27-9  |

## Chemical Structure

