

FMOC-Dap(Boc)-OH *CAS 162558-25-0*

Catalog number: 5110 Unit size: 10 g

Product Details

Storage Conditions Freeze (<-15 °C)

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Solid

Molecular Weight 426.46

Soluble In DMF

Chemical Structure

Applications

FMOC-Dap(Boc)-OH is a common peptide building block. As lysine residue its 3-amino group is often used for carrying a tag moiety such as a fluorescent dye, a quencher or biotin.