

mcKLH *Optimized for Maximum Immunization Response*

Catalog Number: 5607

Unit Size: 10 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	N/A
Soluble In	Water

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Mariculture keyhole limpet hemocyanin (mcKLH) is a purified hemocyanin carrier protein optimized for producing highly effective immunogens with peptide antigens. Due to its high abundance of primary amines, mcKLH facilitates efficient conjugation reactions, while its purified and stabilized form ensures solubility in aqueous solutions. Sourced sustainably from mariculture-grown *Megathura crenulata*, mcKLH exhibits superior immunogenicity attributed to its substantial molecular mass, resulting in a more robust immune response relative to bovine serum albumin (BSA) or ovalbumin. KLH is the carrier protein of choice for the preparation of peptide antigens used in immunization and antibody production, owing to its capacity to elicit a strong immune response when conjugated with small antigens. mcKLH is compatible with hapten-carrier conjugation protocols in both phosphate and MES buffer systems, optimized for NHS-ester and EDC crosslinking methods, respectively. Its substantial structure and abundant reactive groups render mcKLH particularly suited for antibody production and immunological research, ensuring high immunogenicity and minimizing cross-reactivity with mammalian samples.