

Product Information Sheet

Ordering Information

Product Name:	Small Organic Molecule- KLH Conjugate
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving
Molecular Weight:	450-13,000 kD
Chemical Structure:	[Small Organic Molecule] _n – KLH
Soluble in:	Aqueous Buffers

Application Notes

Small Organic Molecule (SOM)–KLH Conjugates are mainly used for developing anti-SOM antibodies that are typically used in an ELISA assay. All our SOM-KLH Conjugates have multiple SOM molecules on each KLH to maximize its immunogenicity. The ratio of SOM/protein is typically kept larger than 15 if SOM does not significantly decrease the water solubility of KLH.