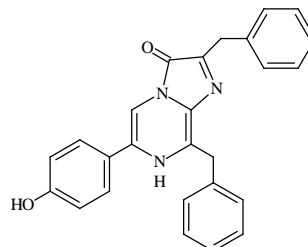


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Aquaphile Coelenterazine *h*, lyophilized solid**CATALOG #:** 10127-50ug/10127**MOLECULAR INFORMATION:** C₂₆H₂₁N₃O₂
Mwt: 407.5**PROPERTIES:**

Color & Form	Yellow solid
Purity	≥ 98% by HPLC
Solubility	Soluble in water and PBS buffer. AVOID DMSO.
Absorption	437 nm
Emission Maxima	466 nm
Extinction Coefficient	9500 (MeOH)

STORAGE AND HANDLING:

Solution is susceptible to oxidation by air. For best results, keep solution from light and store at < -70°C under nitrogen or argon. Keep solid at -20 °C or -70 °C and protect from light under nitrogen or argon for long-term storage. Keep calcium free when stored in solution (avoid using glass container).

APPLICATION:

Aquaphile coelenterazine *h* is a bioluminescent substrate in a special formulation format that is readily dissolved in water or PBS buffer for in vivo use.

TOXICITY:

Unknown

FIRST AID:

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.