



DATA SHEET – Coelenterazine *h*

CHEMILUMINESCENT DETECTION

| Product | Part Number | Quantity | Suggested Reconstitution* |
|-------------------------|-------------|----------|---------------------------|
| Coelenterazine <i>h</i> | 90-0084 | 1 mg | 1 ml Ethanol |
| Coelenterazine <i>h</i> | 90-0084 L | 10 mg | 10 ml Ethanol |

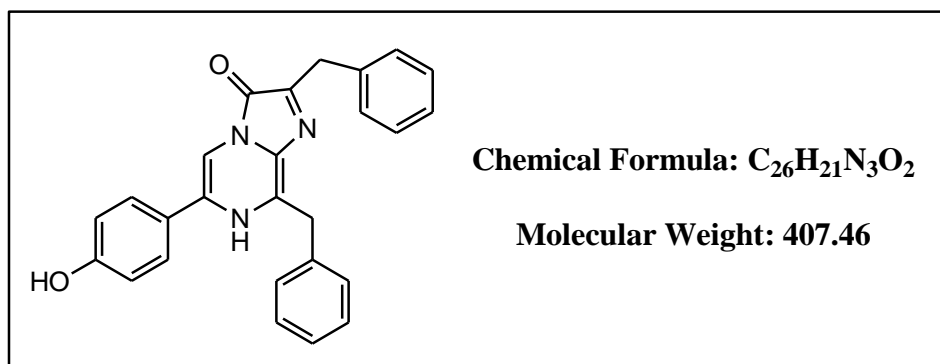
* Reconstitution with the suggested volumes will produce a 2.45 mM solution.

Storage: ≤ -20°C (Protect from light and moisture)

Quality Control Data for Lot Number: 08B0101D

Purity: > 95% by TLC

HPLC: Single peak (using C18RP, CH₃CN/H₂O, at 490nm)



Procedural Notes:

1. To avoid condensation, equilibrate the vial to ambient temperature before opening.
2. Coelenterazine in solution tends to oxidize and gradually decay to Coelenteramide; an oxidation product.
3. Dry Coelenterazine compounds are sparingly soluble in aqueous solutions, and must be dissolved in alcohols, propylene glycol, or straight β-mecaptoethanol prior to making aqueous buffer solutions (Mecaptoethanol may inhibit or destroy luciferase activity without affecting Aqueorin and Obelin; calcium activated photoproteins).
4. Suggested concentration in cells in media is 5uM.

For additional information or Technical Support, contact DiscoverRx or visit the website.

DiscoverRx name and logo are trademarks of DiscoverRx, Corp.

© DiscoverRx Copyrights 2016

Corporate Headquarters

42501 Albrae Street
Fremont, CA 94538 USA
Tel: 510-771-3500 or 866-448-4864 (option 4)
Fax: 510-979-9930
Email: Agreements@discoverx.com
www.discoverx.com

European Regional Headquarters

DiscoverRx Corporation Ltd
Faraday Wharf, Holt Street
Aston Science Park, Birmingham, B7 4BB UK
Phone: +44 121 260 6142
Fax: +44 121 260 6143