

READY-TO-USE COELENTERAZINE

XenoLight RediJect Coelenterazine h

RediJect Coelenterazine h is a bioluminescence substrate in a ready-to-use format developed by the *in vivo* imaging leaders to fit your laboratory workflow. With PerkinElmer's RediJect Coelenterazine you can ensure reproducible results by minimizing variability in formulation, minimizing freeze/thaw and controlling batch to batch variation. RediJect Coelenterazine has been formulated to fit common workflow patterns and optimize results for all your experiments with Renilla luciferase. This ready-to-use substrate formulation allows you to concentrate on animal handling and biology.

- **Pre-formulated** batch controlled Coelenterazine h for *in vivo* use
- **Save time** and effort by minimizing pre-imaging preparation steps
- Dispensed to image **5 animals per vial** (10 vials/kit)
- *In vivo* imaging **quality validated** on IVIS® imaging systems

Part Number: 760506

Molecular Information: C₂₆H₂₁N₃O₂ (MW = 407.5)

Color and Form: Yellow colored solution (Coelenterazine h in Propylene glycol/citrate)

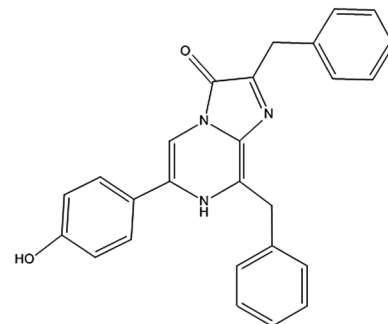
Concentration: 150 µg/mL

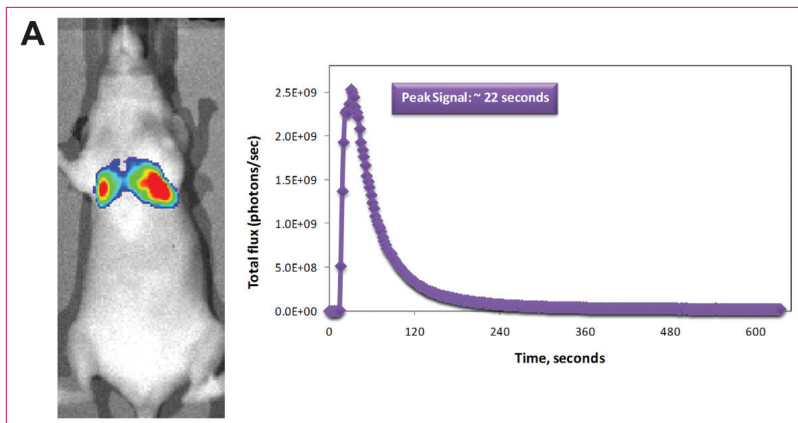
Volume per vial: 10 sterile vials each containing 1.2 mL of 150 µg/mL Coelenterazine h

Storage and Handling: Store at ≤ -20 °C. Just before your experiment, thaw required number of vials in a 37 °C water bath, vortex and it is ready to use. Repeated freeze thaw is not recommended. RediJect Coelenterazine h supports intravenous (i.v.) or intraperitoneal (i.p.) injection.

For intravenous injections, recommended dose is 15 µg/mouse (100 µL/mouse). Load a 1 mL syringe directly from the vial and inject using a 25 gauge needle or higher. Image the mice as soon as possible for peak signal with intravenous injection. For intraperitoneal injection, recommended dose is 30 µg/mouse (200 µL/mouse).

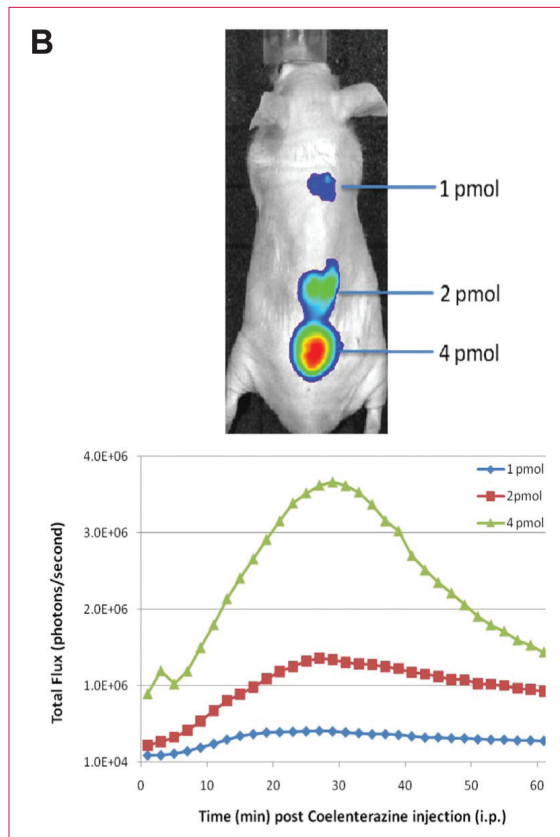
Note: After intravenous injection of RediJect Coelenterazine h mice might experience spasms for a brief period, but should recover shortly.





RediJect Coelenterazine has been optimized for in vivo use with IVIS imaging systems. (A) 5 pmol BRET Qdot 700* were injected intravenously in a nu/nu mouse. 10 minutes post BRET Qdot injection 15 µg/mouse Coelenterazine h was injected intravenously. Chart shows peak signal was observed right after injection at ~ 22 seconds followed by rapid clearance. Images were taken with IVIS Kinetic. (B) 1,2 and 4 pmol BRET Qdot 655 were injected subcutaneously in a nu/nu mouse. 10 minutes post BRET Qdot injection 10 µg/mouse Coelenterazine was injected intraperitoneally. Chart shows that peak signal was reached around 27 minutes post i.p. injection of Coelenterazine.

*BRET Qdot™ 700 were provided by Zymera Inc., 5941 Optical Court, San Jose, CA (<http://zymera.com>) Qdot™ is a registered trademark of Life Technologies, 5791 Van Allen Way, Carlsbad, CA 92008



PerkinElmer *in vivo* imaging reagents are intended for animal research and not for use in humans.

Learn more at www.perkinelmer.com/invivoreagents

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

010429_01

Printed in USA