



**EAGLE**  
BIOSCIENCES

**Glucuronidase Sample Treatment Kit**

Catalog Number: ISO34-PREP

Size: 2 vials, lyophilized

Storage: -20°C

**PRODUCT SPECIFICATION SHEET**

- DESCRIPTION:** Isoprostanes are primarily generated by free radical mechanisms making them excellent biomarkers for systemic oxidant stress. It is now known that isoprostanes are extensively metabolized in humans. Eagle Biosciences has found that an average of 50% of the isoprostane excreted in human urine is conjugated to glucuronic acid. The extent of glucuronidation among individuals ranges significantly from 28% to 80%. In light of this information it is strongly recommended that specimens be pretreated with  $\beta$ -glucuronidase prior to isoprostane analysis to provide a more accurate assessment of oxidative stress.
- This Glucuronidase Sample Treatment Kit provides sufficient materials and methods for the treatment of 40 samples allowing the user to differentiate the inter-individual differences in glucuronidation and measure the total systemic isoprostane output.
- CONCENTRATION:** 125,000 U/mL/vial upon reconstitution
- PRESENTATION:** 22,500 U/vial lyophilized powder
- STORAGE:** -20°C. This product is intended for use in its entirety once reconstituted - **AVOID REPEATED FREEZE-THAW CYCLES.**
- APPLICATIONS:** For use in the removal of conjugated glucuronide from isoprostane.
- SPECIES SPECIFICITY:** Not species specific.
- DIRECTIONS FOR USE:**
- 1) Reconstitute powder with 180  $\mu$ L of deionized water and place on ice for no longer than one hour prior to use. Due to the viscosity of the reconstituted solution, take care to avoid bubbles.
  - 2) For every 100  $\mu$ L of stock urine to be assayed, add 8  $\mu$ L of Glucuronidase. Seal container and mix solution by inversion.
  - 3) Incubate the mixture at 37°C for 2 hours.
  - 4) Your sample is now ready for dilution and assay.

*Product Developed and Manufactured in the USA*

---

EAGLE BIOSCIENCES, INC.  
20A NW Blvd., Suite 112  
Nashua, NH 03063

Phone: 617-419-2019  
Fax: 617-419-1110  
Email: [info@EagleBio.com](mailto:info@EagleBio.com)  
[www.EagleBio.com](http://www.EagleBio.com)