

3D InSight™ Krebs-Ringer HEPES Buffer

3D InSight™ Krebs-Ringer HEPES Buffer is a physiological buffer designed for the assessment of hormone secretion from 3D InSight™ Islet Microtissues.

Key Features

- Formulated for optimal evaluation of pancreatic β -cell and α -cell function
- Compatible with commercially available ELISAs for insulin, proinsulin and glucagon
- Maintains ideal physiological pH range, intra- and extracellular osmotic balance during assays.

3D InSight™ Krebs-Ringer HEPES Buffer

Part Number: CS-07-051-01

- Certified for the assessment of
 - Insulin Secretion (TP013)
 - Glucagon Secretion (TP015)
- Available only when purchased with 3D InSight™ Islet Microtissues



Preparation Instructions

1. Prepare KRHB working solution by adding 0.5% (w/v) Bovine Serum Albumin (BSA, RIA grade, low insulin, recommended: A7888 Sigma-Aldrich).
2. Gently shake until BSA is completely dissolved.
3. Measure pH, which should remain at 7.4 after the addition of BSA.
4. Filter the solution through a 0.2 μ m bottle top filter.
5. Store at 4° C for up to 2 months after the addition of BSA.
6. Freshly add the required amount of glucose prior to secretion experiments.

Related Products

- 3D InSight™ Islet Microtissues (human and rat)
- 3D InSight™ Islet Maintenance Medium
- 3D InSight™ Islet Assay Medium
- ALPCO STELLUX® Chemi Insulin ELISAs

IS-500-032-01-A

InSphero AG
Schlieren, Switzerland
© +41 44 515 04 90

InSphero Inc.
Brunswick, ME, USA
© +1 800-779-7558

Email and Website
info@insphero.com
www.insphero.com

InSphero is ISO 9001:2015 certified
All rights reserved, © 2018 InSphero AG. 3D InSight and Akura are trademarks of InSphero AG. All other trademarks cited are the property of their respective owners. For life science research only. Not for use in diagnostic procedures.