

HTRF[®] Insulin mouse serum kit



THE SENSITIVITY, DYNAMIC RANGE, AND RELIABILITY OF ELISA. A FRACTION OF THE TIME AND SAMPLE VOLUME.

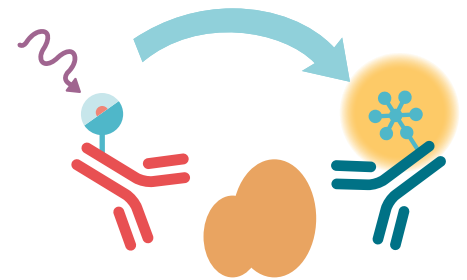
For many years, researchers have sought the ideal insulin assay combining sensitivity, accuracy, low sample consumption, and ease-of-use. Their search has now ended.

The Cisbio mouse serum insulin assay offers superior sensitivity and dynamic range to quantify insulin reliably in mouse serum and plasma samples. Validated on metabolic and diabetic mouse research models, this kit offers outstanding correlation with ELISA assays.

Built on a truly homogeneous protocol, the assay does not require any washing or complex extraction steps, which saves significant time as compared to ELISA methods.

Designed for *in vivo* experiments on mouse models secreting all insulin levels.

- High sensitivity
- Broad dynamic range
- Thoroughly validated
- Highly correlated with ELISA
- Protocol available in 5 μ L or 2 μ L samples
- Truly homogeneous time-saving technology



Insulin is detected in a sandwich assay using two different specific monoclonal antibodies: one labeled with cryptate (donor) and the second with XL665 (acceptor).

Reagents are lyophilized for increased stability

DIABETES

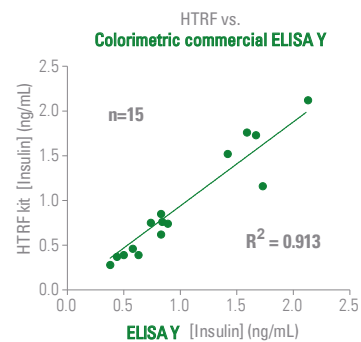
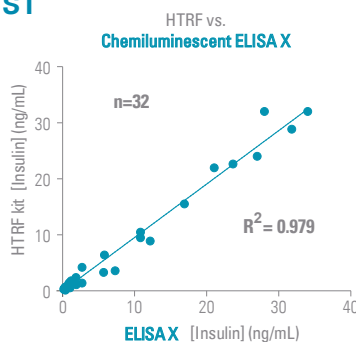
ASSAY CHARACTERISTICS: OUR SENSITIVITY SERVED BY PRECISION AND DYNAMIC RANGE

Sample volume	2 or 5 µL
Assay range	0.24 - 8 ng/mL
The LoD detection limit is calculated as two standard deviations above the standard 0. The HTRF® Insulin assay is calibrated against the NIBSC international standard 66/304. 1ng Insulin HTRF® ↔ 1ng Insulin NIBSC 66/304 (i.e.: 2.3x10 ⁻⁰⁵ IU).	0.064 ng/mL (LoD obtained for 5 µL sample)
Incubation time	Overnight at RT
Recovery upon addition	95 - 117 % (104 %)
Recovery upon dilution	98 - 118 % (109 %)
Intra-assay (n=24) on 5 mouse serum samples	4 - 7.6% (5.1%)
Inter-assay (n=4) on 5 mouse serum samples	4 - 14.6% (9.1%)

HIGH CORRELATION FOR RESULTS YOU CAN TRUST

A study of 32 and 15 normal, diabetic, and obese mouse plasma samples using two commercially available ELISA kits shows strong correlation with the HTRF insulin assay.

Results shows the insulin quantification obtained in each individual assay.



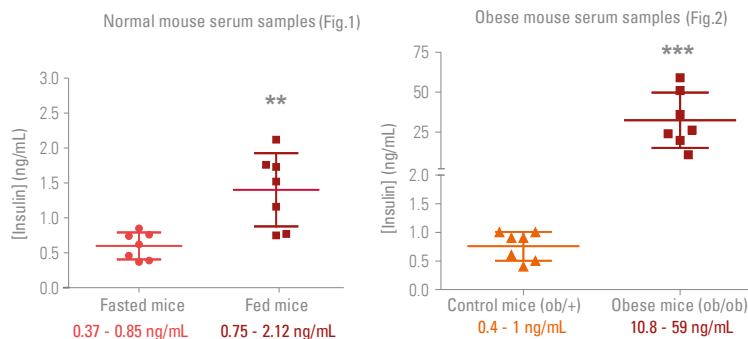
THOROUGHLY VALIDATED ON COMMONLY USED METABOLIC MOUSE MODELS

Fig.1 shows Insulin levels in C57BL/6J mice randomized into two groups of 7 animals. One group was fasted overnight, while the second was fed normally.

Fig.2 shows Insulin levels in one group of eight ob/ob obese mice and one group of eight ob/+ control mice.

Mice were sacrificed and the blood was collected through cardiac puncture for serum preparation.

Results shows the insulin quantification obtained in each individual assay using the HTRF Mouse serum Insulin kit reagents.



Samples were kindly provided by JF Tanti's research team: Cellular and Molecular Physiopathology of Obesity and Diabetes, UMR INSERM U1065/UNS, C3M, Nice, France

ORDERING INFORMATION

PRODUCT	PART#
Insulin Mouse Serum kit - 200 tests*	62IN3PEF
Insulin Mouse Serum kit - 5 x 200 tests*	62IN3PEB

* based on 20 µL assay volume

GET EXACTLY WHAT YOU NEED WITH OUR FULL SERVICE CAPABILITIES



Discover a partner dedicated to meeting your exact needs, from our assay chemistry to the collaboration-oriented nature of our relationship with customers. Cisbio is firmly committed to the belief that "interaction is everything." Talk to us and get the right tools for your work.

Keep your diabetes study moving forward with faster, more robust, and physiologically relevant results.
Call us or visit www.htrf.com

FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567

China +86 10 8080 9288 Japan +81-(0)43-306-8712

Visit www.cisbio.com to find a list of our regional distributors



cisbio