

Biomarker and Cell Signaling Assays for Fibrosis and NASH



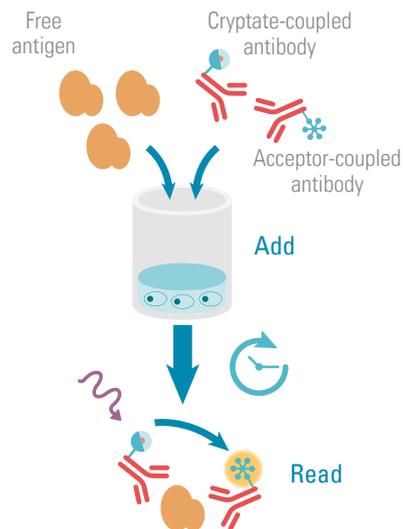
HTRF® SOLUTIONS FOR NASH

Nonalcoholic steatohepatitis (NASH) is a specific type of nonalcoholic fatty liver disease (NAFLD) where fat accumulation, or steatosis, occurs in the liver due to non-alcohol related causes. While the precise mechanisms of the disease are unknown, steatosis has been linked to insulin resistance and can result in the stimulation of hepatic stellate cells, leading to fibrosis, inflammation, and eventually cirrhosis.

Cisbio offers advanced solutions for evaluating NASH mechanisms and treatments with HTRF assays.

Elucidate NASH mechanisms with Cisbio:

- Robust and high-performing assay portfolio for the investigation of NASH molecular mechanisms
- Specific cell type & multiple cell type products
- Dedicated scientific consultants provide support and experiment troubleshooting



HTRF Technology, a powerful, easy alternative to ELISA & Western Blot

- Easy add-and-read protocol saves you time with NO washing steps
- Low sample volume consumption saves laboratory resources

HEPATOCTE CELLS

| PRODUCT | PART# |
|------------------------------------|-------------|
| ACC phospho-S79 kit | 64ACCPEG |
| AMPK phospho -T172 * | 64MPKPEG |
| BAD phospho-S112 kit * | 64BADPEG |
| CREB phospho-S133 kit * | 63ADK052PEG |
| EIF2 alpha phospho-S52 kit * | 64EF2PEG |
| IRS1 phospho-S307(m)/S312(h) kit * | 64IRSPEG |
| IRS2 total kit | 64IRTPEG |
| PARP cleaved | 64PARPEG |

IMMUNE CELLS

| PRODUCT | PART# |
|-----------------------|-------------|
| cAMP-Gi | 62AM9PEG |
| cRAF phospho-S338 kit | 63ADK018PEG |
| IP-One Gq | 62IPAPEB |

HEPATIC STELLATE CELLS (HSCs)

| PRODUCT | PART# |
|---------------------------------------|-------------|
| Alpha-SMA | 62ASMAPEG |
| Beta Catenin phospho-T41/S33/37 kit * | 64CATPEG |
| Hyaluronic Acid | 63ADK089PEG |
| MMP2 human kit | 62MM9PEG |
| MMP9 human kit | 62MM9PEG |
| P3NP ELISA | P3NP-EL-US |
| PDGFR-beta phospho-Y751 kit | 63ADK040PEG |
| Pro-Collagen Type 1 | 63ADK014PEG |
| TIMP1 human kit | 64TIMPEG |
| YAP phospho-S127 kit * | 64YAPPEG |

Part numbers reflect the smallest available assay format.

* Also available in total kit

PRO-INFLAMMATORY CYTOKINES & CHEMOKINES

| PRODUCT | PART# |
|----------------------|-------------|
| CCL2 human | 62HCCL2PEG |
| CCL5 human kit | 62HCCL5PEG |
| IFN alpha human kit | 62HIFNAPEG |
| IFN beta human kit | 62HIFNBPEG |
| IL1 beta human | 62HIL1BPET |
| IL1 beta mouse | 62MIL1PEG |
| IL12 human | 62HIL12PEG |
| IL17 human | 62HIL17PEG |
| IL17 mouse | 62MIL17PEG |
| IL18 human | 62HIL18PEG |
| IL6 human | 62HIL06PET |
| IL6 mouse | 62MIL06PEG |
| IL8 human kit | 62HIL08PEG |
| TGF beta | 62HTGBPEG |
| TNF alpha cynomolgus | 62CTNFAPEG |
| TNF alpha human | 62HTNFAPET |
| TNF alpha mouse | 62MNTNFAPEG |

APPLICABLE TO MULTIPLE CELL TYPES

| PRODUCT | PART# |
|----------------------------------|-------------|
| AKT phospho -S473 kit * | 64AKTPEG |
| ATF2 phospho-T711 kit | 63ADK015PEG |
| cJun phospho-S63 kit | 64JUNPEG |
| ERK phospho-T202/Y204 kit * | 64AERPEG |
| FAK phospho-Y397 kit * | 64FAKPEG |
| FOXO1 phospho -S256 kit * | 64FOXPEG |
| GSK3 alpha phospho-S21 kit * | 64GPAPEG |
| GSK3 beta phospho-S9 kit * | 64GPBPEG |
| HIF-1alpha kit | 64HIFPEG |
| IRF3 phospho-S386 kit * | 6FRF3PEG |
| JNK phospho-T183/Y185 kit | 64JNKPEG |
| MEK1 phospho S218/22 kit * | 64ME1PEG |
| MEK1 phospho-S298 kit * | 63ADK201PEG |
| MKK4 phospho-S257 kit NEW | 64MK4PEG |
| mTOR phospho-S2448 kit | 64TORPEG |
| NFkB phospho-S536 kit | 64NFBPEG |
| P38 phospho-T180/Y182 kit | 64P38PEG |
| SMAD2 phospho-S465/467 kit * | 63ADK102PEG |
| SMAD3 phospho-S423/425 kit * | 63ADK025PEG |
| STAT3 phospho-Y705 kit | 62AT3PEG |
| TBK1 phospho-S172 kit | 64TBKPEG |

EXPERT SERVICES TO MEET YOUR PRECISE NEEDS

In addition to providing ready-to-use kits with all the necessary reagents, Cisbio's team of experienced experts works hand-in-hand with our clients to create custom solutions. When you have specific requirements for fibrosis and NASH, Cisbio is the ideal trusted partner to develop the assays you need.

Keep your therapeutic research moving forward with faster, more robust, physiologically relevant results. Call us or visit www.cisbio.com



FOR MORE INFORMATION

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

China +86 21 5018 9880 Japan +81-(0)43-306-8712

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



cisbio