

Biomarker and Cell Signaling Assays for Fibrosis and NASH



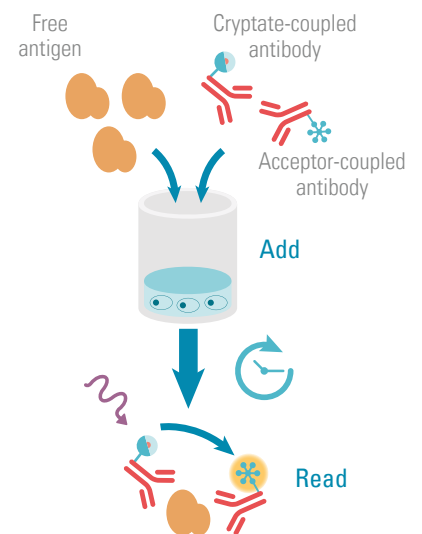
HTRF® SOLUTIONS FOR NASH

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	62IPAPEG

HEPATIC STELLATE CELLS (HSCs)

PRODUCT	PART#
Alpha-SMA	62ASMAPEG
Beta Catenin phospho-T41/S33/37 kit *	64CATPEG
Hyaluronic Acid	63ADK089PEG
Hyaluronic Acid ELISA	HYAL-US
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
mTOR phospho-S2448 kit	64TORPEG
NFkB phospho-S536 kit	64NFBPEG
P38 phospho-T180/Y182 kit	64P38PEG
SMAD2 phospho-S465/467 kit * NEW	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

