



LANCE *Ultra*

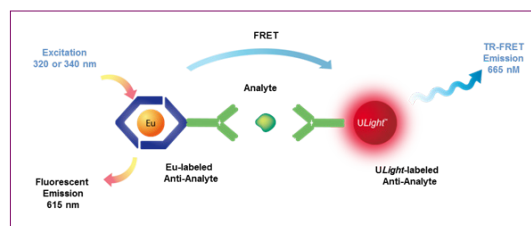
TR-FRET Biomarker
Detection Kits

The original, proven TR-FRET technology that can brighten today and free up tomorrow

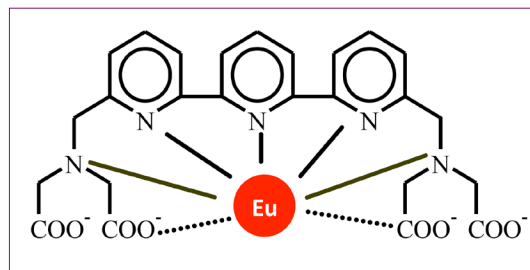
From assay development to primary and secondary screening, LANCE® *Ultra* Technology provides researchers with robust and easy-to-use assays adaptable to a variety of targets on a broad range of multimode detectors, with the fastest time-to-results in a commercially-available TR-FRET assay.

- No-wash/separation steps
- Homogeneous dispense-and-read technology
- Sensitive detection
- High reproducibility
- Faster time-to-results
- Easy automation
- 384-well and 1536-well formats

LANCE and LANCE (LANThanide Chelate Excitation) *Ultra* are our TR-FRET (time-resolved fluorescence resonance energy transfer), homogeneous (no wash) technologies. One binding partner of interest is labeled with a donor fluorophore (a LANCE Europium chelate) and a second one is labeled with an acceptor fluorophore (*ULight*™ dye). Upon excitation at 320–340 nm, energy can be transferred from the donor Europium chelate to the acceptor fluorophore if sufficiently close for FRET (~10 nm). This results in the emission of light detectable at 665 nm.

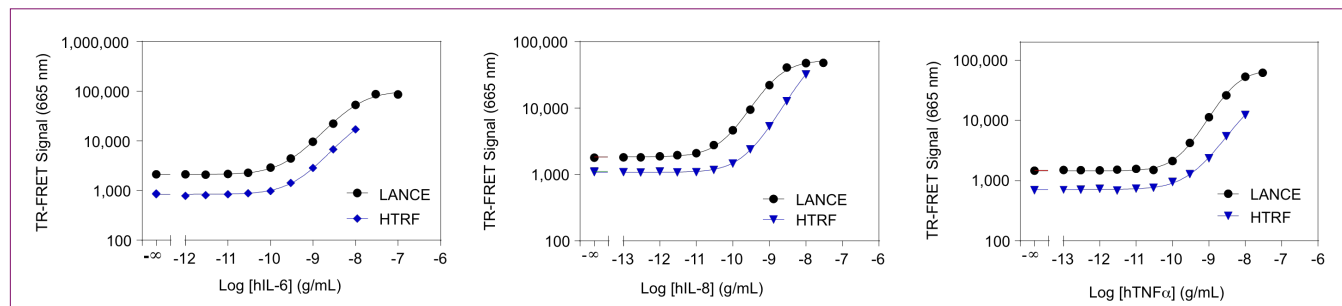


One antibody of interest is labeled with a donor fluorophore (a LANCE Europium chelate) and the second molecule is labeled with an acceptor fluorophore (*ULight*™ dye). Upon excitation at 320 or 340 nm, energy can be transferred from the donor Europium chelate to the acceptor fluorophore if sufficiently close for FRET (~10 nm). This results in the emission of light at 665 nm.



Representative LANCE chelate structure. Stable, hydrophilic, and no potassium fluoride (KF) needed.

Comparison with HTRF Assays



Comparison of LANCE *Ultra* and HTRF® Assays. Standard curves obtained for IL-6, IL-8, and TNFα after one-hour incubation. Sensitivity for IL-6 (LDL=10-15 pg/ml) and TNFα (LDL=20 pg/ml) kits were equivalent for LANCE *Ultra* and HTRF® products, whereas LANCE *Ultra* IL-8 (LDL=5 pg/ml) kits showed slightly superior sensitivity for detecting IL-8 compared to HTRF® (LDL=20 pg/ml) kits.

Latest LANCE *Ultra* Biomarker Detection Kits

NAME OF KIT	500 ASSAY POINTS	10,000 ASSAY POINTS
ANGIOGENESIS		
LANCE <i>Ultra</i> ANGPT1 Kit	TRF3014C	TRF3014M
LANCE <i>Ultra</i> ANGPTL3 Kit	TRF3016C	TRF3016M
LANCE <i>Ultra</i> ANGPTL4 Kit	TRF3017C	TRF3017M
ANIMAL HEALTH		
LANCE <i>Ultra</i> IgG (bovine) Kit	TRF1530C	TRF1530M
AUTOPHAGY		
LANCE <i>Ultra</i> LC3B Kit	TRF1306C	TRF1306M
BIOLOGICS / BIOPROCESS		
LANCE <i>Ultra</i> human IgA Kit	TRF1262C	TRF1262M
LANCE <i>Ultra</i> human IgE Kit	TRF1292C	TRF1292M
LANCE <i>Ultra</i> human IgG Kit	TRF1205C	TRF1205M
LANCE <i>Ultra</i> human IgM Kit	TRF1263C	TRF1263M
LANCE <i>Ultra</i> mouse IgG Kit	TRF1529C	TRF1529M
LANCE <i>Ultra</i> rabbit IgG Kit	TRF1104C	TRF1104M
LANCE <i>Ultra</i> human Kappa Light Chain Kit	TRF3023C	TRF3023M
LANCE <i>Ultra</i> human Lambda Light Chain Kit	TRF3024C	TRF3024M
LANCE <i>Ultra</i> c-Myc Kit	TRF1111C	TRF1111M
LANCE <i>Ultra</i> CHO HCP Kit	TRF1301C	TRF1301M
LANCE <i>Ultra</i> CHO-P Kit	TRF1210C	TRF1210M
LANCE <i>Ultra</i> E.coli HCP Kit	TRF1261C	TRF1261M
LANCE <i>Ultra</i> PER.C6 HCP Kit	TRF1302C	TRF1302M
LANCE <i>Ultra</i> Residual DNA Kit	TRF1331C	TRF1331M
LANCE <i>Ultra</i> Residual Protein A Kit	TRF1287C	TRF1287M
CANCER		
LANCE <i>Ultra</i> Bcl-2 Kit	TRF1384C	TRF1384M
LANCE <i>Ultra</i> COL1A1 Kit	TRF1371C	TRF1371M
LANCE <i>Ultra</i> COX-2 Kit	TRF1376C	TRF1376M
LANCE <i>Ultra</i> E-cadherin Kit	TRF1370C	TRF1370M
LANCE <i>Ultra</i> EGF Kit	TRF1339C	TRF1339M
LANCE <i>Ultra</i> EGFR Kit	TRF1340C	TRF1340M
LANCE <i>Ultra</i> ErbB2 Kit	TRF1235C	TRF1235M
LANCE <i>Ultra</i> Fibronectin Kit	TRF1351C	TRF1351M
LANCE <i>Ultra</i> MANF Kit	TRF3009C	TRF3009M
LANCE <i>Ultra</i> MMP-3 Kit	TRF1284C	TRF1284M
LANCE <i>Ultra</i> MMP-7 Kit	TRF1378C	TRF1378M
LANCE <i>Ultra</i> MMP-8 Kit	TRF1374C	TRF1374M
LANCE <i>Ultra</i> MMP-12 Kit	TRF1375C	TRF1375M
LANCE <i>Ultra</i> Nidogen-1 Kit	TRF1367C	TRF1367M
LANCE <i>Ultra</i> Oncostatin M Kit	TRF1341C	TRF1341M
LANCE <i>Ultra</i> PD-1 Kit	TRF1343C	TRF1343M
LANCE <i>Ultra</i> PD-L1 Kit	TRF1355C	TRF1355M
LANCE <i>Ultra</i> Perlecan Kit	TRF1377C	TRF1377M
LANCE <i>Ultra</i> PIP Collagen Kit	TRF1353C	TRF1353M
LANCE <i>Ultra</i> PTEN Kit	TRF1380C	TRF1380M
LANCE <i>Ultra</i> VEGF (mouse) Kit	TRF1520C	TRF1520M
LANCE <i>Ultra</i> VEGF Kit	TRF1201C	TRF1201M
CARDIOVASCULAR		
LANCE <i>Ultra</i> Fibrinogen Kit	TRF1352C	TRF1352M
LANCE <i>Ultra</i> tPA Kit	TRF1250C	TRF1250M
LANCE <i>Ultra</i> Transferrin Kit	TRF1311C	TRF1311M
CENTRAL NERVOUS SYSTEM		
LANCE <i>Ultra</i> A β (1-40) Kit	TRF1345C	TRF1345M
LANCE <i>Ultra</i> ApoE Kit	TRF1395C	TRF1395M
LANCE <i>Ultra</i> BDNF Kit	TRF1317C	TRF1317M
LANCE <i>Ultra</i> Frataxin Kit	TRF1322C	TRF1322M

NAME OF KIT	500 ASSAY POINTS	10,000 ASSAY POINTS
LANCE <i>Ultra</i> GAD67 Kit	TRF1397C	TRF1397M
LANCE <i>Ultra</i> Tau Kit	TRF1271C	TRF1271M
LANCE <i>Ultra</i> TDP-43 Kit	TRF1387C	TRF1387M
INFLAMMATION/CYTOKINE		
LANCE <i>Ultra</i> CCL2/MCP-1 Kit	TRF1244C	TRF1244M
LANCE <i>Ultra</i> CCL2/MCP-1 α (mouse) Kit	TRF1509C	TRF1509M
LANCE <i>Ultra</i> CCL3/MIP1 α Kit	TRF1257C	TRF1257M
LANCE <i>Ultra</i> CCL4 Kit	TRF1258C	TRF1258M
LANCE <i>Ultra</i> CCL5 Kit	TRF1289C	TRF1289M
LANCE <i>Ultra</i> CRP Kit	TRF1233C	TRF1233M
LANCE <i>Ultra</i> CXCL1/GRO- α Kit	TRF1279C	TRF1279M
LANCE <i>Ultra</i> CXCL9 Kit	TRF1280C	TRF1280M
LANCE <i>Ultra</i> CXCL10 Kit	TRF1259C	TRF1259M
LANCE <i>Ultra</i> IFN α Kit	TRF1297C	TRF1297M
LANCE <i>Ultra</i> IFN γ Kit	TRF1217C	TRF1217M
LANCE <i>Ultra</i> IFN γ (mouse) Kit	TRF1501C	TRF1501M
LANCE <i>Ultra</i> IFN- λ 2 Kit	TRF3039C	TRF3039M
LANCE <i>Ultra</i> IL-1 α Kit	TRF1238C	TRF1238M
LANCE <i>Ultra</i> IL-1 β Kit	TRF1220C	TRF1220M
LANCE <i>Ultra</i> IL-1RA Kit	TRF1358C	TRF1358M
LANCE <i>Ultra</i> IL-1 β (mouse) Kit	TRF1503C	TRF1503M
LANCE <i>Ultra</i> IL-2 Kit	TRF1221C	TRF1221M
LANCE <i>Ultra</i> IL-4 Kit	TRF1266C	TRF1266M
LANCE <i>Ultra</i> IL-5 Kit	TRF1267C	TRF1267M
LANCE <i>Ultra</i> IL-6 Kit	TRF1223C	TRF1223M
LANCE <i>Ultra</i> IL-6 (mouse) Kit	TRF1504C	TRF1504M
LANCE <i>Ultra</i> IL-8 Kit	TRF1224C	TRF1224M
LANCE <i>Ultra</i> IL-11 Kit	TRF1293C	TRF1293M
LANCE <i>Ultra</i> IL-12/p40 Kit	TRF1382C	TRF1382M
LANCE <i>Ultra</i> IL-17A Kit	TRF1219C	TRF1219M
LANCE <i>Ultra</i> IL-17A (mouse) Kit	TRF1514C	TRF1514M
LANCE <i>Ultra</i> IL-17F Kit	TRF1391C	TRF1391M
LANCE <i>Ultra</i> IL-23A Kit	TRF3006C	TRF3006M
LANCE <i>Ultra</i> IL-28B Kit	TRF1365C	TRF1365M
LANCE <i>Ultra</i> IL-31 Kit	TRF3003C	TRF3003M
LANCE <i>Ultra</i> TNF α Kit	TRF1208C	TRF1208M
LANCE <i>Ultra</i> TNF α (mouse) Kit	TRF1505C	TRF1505M
METABOLIC		
LANCE <i>Ultra</i> Adiponectin Kit	TRF1209C	TRF1209M
LANCE <i>Ultra</i> Albumin Kit	TRF1294C	TRF1294M
LANCE <i>Ultra</i> ApoA1 Kit	TRF1389C	TRF1389M
LANCE <i>Ultra</i> Apo-CIII Kit	TRF1396C	TRF1396M
LANCE <i>Ultra</i> C-peptide (mouse/rat) Kit	TRF1510C	TRF1510M
LANCE <i>Ultra</i> Chorionic Gonadotropin Kit	TRF3001C	TRF3001M
LANCE <i>Ultra</i> Cortisol Kit	TRF1314C	TRF1314M
LANCE <i>Ultra</i> FSH Kit	TRF1394C	TRF1394M
LANCE <i>Ultra</i> GLP-1 (7-36) Kit	TRF1215C	TRF1215M
LANCE <i>Ultra</i> Glucagon Kit	TRF1312C	TRF1312M
LANCE <i>Ultra</i> IGF-1 Kit	TRF1236C	TRF1236M
LANCE <i>Ultra</i> IGF-1 (rat) Kit	TRF1552C	TRF1552M
LANCE <i>Ultra</i> Insulin Kit	TRF1204C	TRF1204M
LANCE <i>Ultra</i> Leptin Kit	TRF1225C	TRF1225M
LANCE <i>Ultra</i> Leptin (mouse) Kit	TRF1521C	TRF1521M
LANCE <i>Ultra</i> MBNL1 Kit	TRF3008C	TRF3008M
LANCE <i>Ultra</i> PAI-1 Kit	TRF1286C	TRF1286M
LANCE <i>Ultra</i> PCSK9 Kit	TRF1270C	TRF1270M
LANCE <i>Ultra</i> TSH Kit	TRF3005C	TRF3005M

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2016-2017 PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.