

LANCE *Ultra* JMJD3 Histone H3-Lysine 27 Demethylase Assay

U-TRF #42

LANCE® *Ultra*

Authors

Anne Labonté
Mathieu Arcand
Julie Blouin
Mireille Caron
Claire Normand
Lucille Beaudet
Jaime Padrós

PerkinElmer, Inc.
Montreal, QC
Canada, H3J 1R4

This LANCE *Ultra* immunodetection assay measures the demethylation of a biotinylated Histone H3 (21-44) peptide tri-methylated at lysine 27.

Europium-anti-di/mono-methyl-Histone H3 Lysine 27 (H3K27me2-1) Antibody

- TRF0406-D: 10 µg, 1,562 assay points*
- TRF0406-M: 100 µg, 15,625 assay points*

*40 fmol/assay point

Peptidic Substrate Sequence:

ATKAARK(me3)SAPATGGVKKPHRYRP-GG-K(Biotin)-OH

LANCE *Ultra* Assays

LANCE *Ultra* time-resolved fluorescence resonance energy transfer (TR-FRET) assays use a proprietary europium chelate donor dye, W1024 (Eu), together with *ULight*™, a small molecular weight acceptor dye with a red-shifted fluorescent emission.

In this technical note, we present the optimization of an epigenetic enzymatic assay using a biotinylated Histone H3-derived peptide as substrate. The modified peptide is captured by the Eu-labeled antibody (Eu-Ab) and *ULight*-Streptavidin (*ULight*-SA), which brings the Eu donor and *ULight* acceptor dye molecules into close proximity. Upon irradiation at 320 or 340 nm, the energy from the Eu donor is transferred to the *ULight* acceptor dye which, in turn, generates light at 665 nm. The intensity of the light emission is proportional to the level of biotinylated substrate modification.

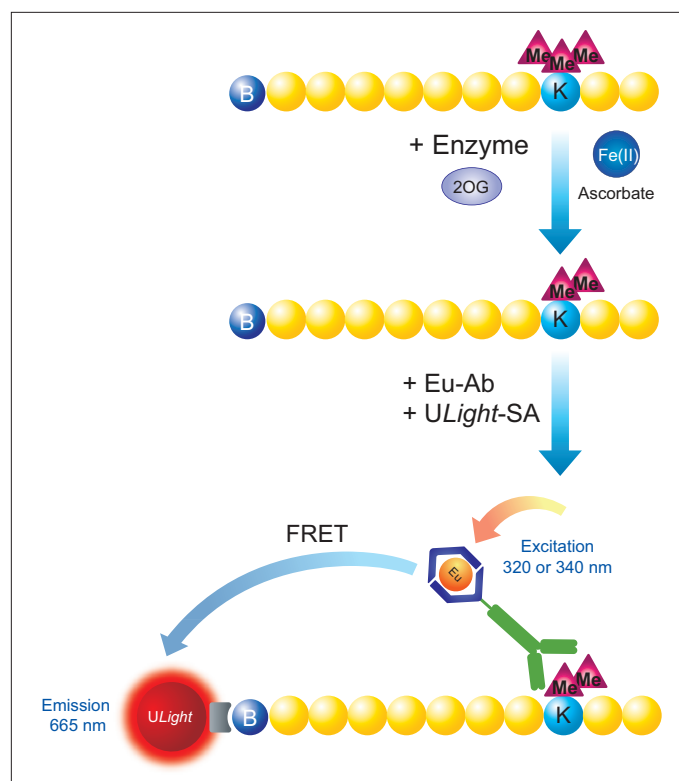


Figure 1. Schematic representation of the LANCE *Ultra* detection of a modified histone peptide.

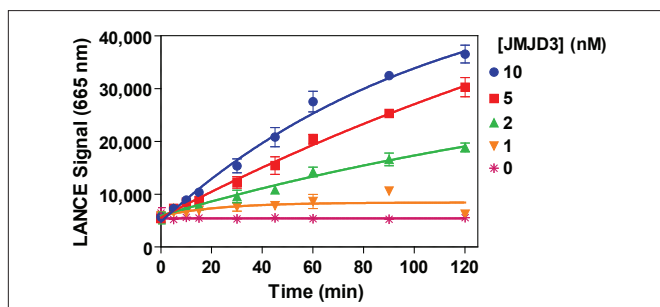
Development of a JMJD3 Histone H3-Lysine 27 Demethylase Assay:

Reagents needed for the assay:

Europium-anti-di/mono-methyl-Histone H3 Lysine 27 (H3K27me2-1) Antibody	PerkinElmer # TRF0406
LANCE Ultra ULight-Streptavidin	PerkinElmer # TRF0102
Histone H3 (21-44), H3K27(me3) peptide, biotinylated	AnaSpec # 64367
LANCE Detection Buffer, 10X	PerkinElmer # CR97-100
JMJD3 (human), recombinant	BPS BioScience # 50115
White opaque OptiPlate™-384	PerkinElmer # 6007299
TopSeal™-A films	PerkinElmer # 6005185
α-Ketoglutaric acid potassium salt (2OG)	Sigma # K2000
(+) Sodium L-ascorbate	Sigma # 11140
Ammonium iron(II) sulfate hexahydrate (Fe(II))	Sigma # 215406
2,4-Pyridinedicarboxylic acid (2,4-PDCA)	Sigma # P63395
Ethylenediaminetetraacetic acid (EDTA)	Invitrogen #15575-038

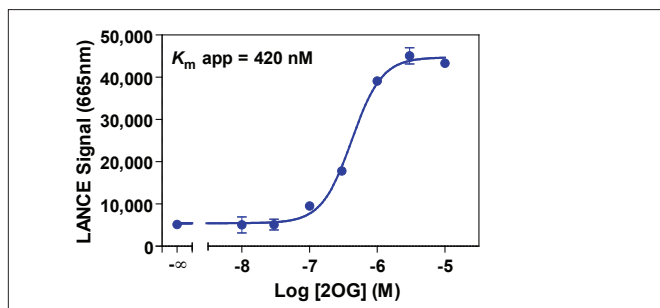
Assay Buffer: 50 mM HEPES pH 7.5, 0.01% Tween-20 and 0.01% BSA

Experiment 1: Enzyme Titration and Time-Course



Enzymatic progress curves were performed by incubating JMJD3 at concentrations ranging from 1 to 10 nM with 200 nM biotinylated Histone H3K27me3 peptide substrate plus 50 μM 2OG, 5 μM Fe(II) and 100 μM ascorbate. Reactions were stopped by the addition of EDTA at indicated times. The Detection Mix was then added and signal was read after 60 min. A 120 min reaction time using 5 nM enzyme was selected for all subsequent experiments.

Experiment 2: 2OG Titration



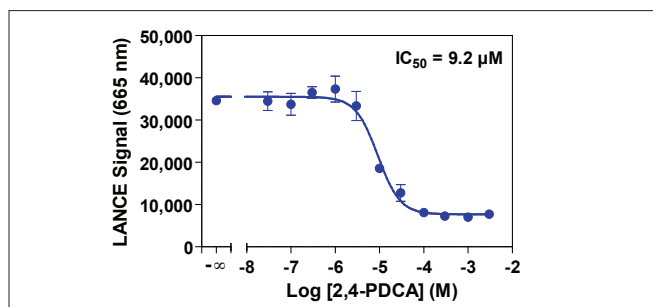
Serial dilutions of 2OG ranging from 10 nM to 10 μM were added to 5 nM JMJD3 and 200 nM biotinylated Histone H3K27me3 peptide substrate plus 5 μM Fe(II) and 100 μM ascorbate. A 500 nM 2OG concentration was selected for subsequent experiments.

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

Standard Protocol

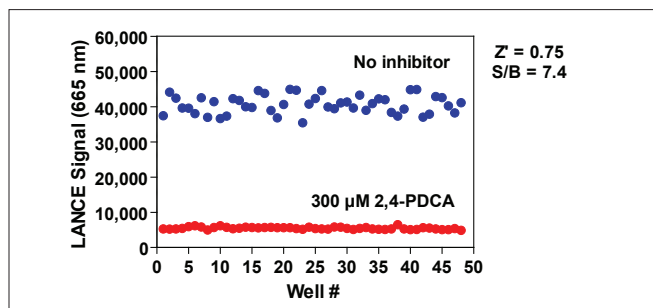
- Dilute JMJD3 enzyme, 2OG, Fe(II), ascorbate, 2,4-PDCA (inhibitor) and biotinylated peptide substrate in Assay Buffer just before use.
- Add to the wells of a white OptiPlate-384:
 - 5 μL of inhibitor (2X) or Assay Buffer
 - 2.5 μL of enzyme (4X)
 - 2.5 μL of biotinylated Histone H3K27me3 peptide/2OG/Fe(II)/ascorbate mix (4X).
 For 2OG titration, add 2OG dilutions independently of substrate.
- Cover the plate with TopSeal-A film and incubate at room temperature (RT).
- Prepare a 4X Detection Mix by diluting the Eu-Ab to 8 nM and ULight-Streptavidin to 200 nM in 1X LANCE Detection Buffer (final concentrations of 2 nM and 50 nM, respectively, in 20 μL total assay volume).
- Prepare a 4X Stop Solution containing 4 mM EDTA in 1X LANCE Detection Buffer (final concentration of 1 mM in 20 μL total assay volume).
 - 5 μL of EDTA Stop Solution and incubate 5 min at RT
 - 5 μL of Detection Mix
- Cover with TopSeal-A film and incubate for 60 min at RT.
- Remove the TopSeal-A film and read signal with the EnVision® Multilabel Reader in TR-FRET mode (excitation at 320 or 340 nm & emission at 665 nm).

Experiment 3: Enzyme Inhibition



Serial dilutions of 2,4-PDCA ranging from 30 nM to 3 mM were pre-incubated for 15 min with 5 nM JMJD3. Enzymatic reactions were initiated by the addition of 200 nM biotinylated Histone H3K27me3 peptide substrate plus 500 nM 2OG, 5 μM Fe(II) and 100 μM ascorbate. Enzymatic reactions contain 1% DMSO.

Experiment 4: Z'-factor Determination



JMJD3 (5 nM) was incubated with or without 300 μM 2,4-PDCA. Enzymatic reactions were initiated by the addition of 200 nM biotinylated Histone H3K27me3 peptide substrate plus 500 nM 2OG, 5 μM Fe(II) and 100 μM ascorbate. Enzymatic reactions contain 0.05% BSA and 1% DMSO.



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

Copyright © 2010-2011, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.