

DNMT Colorimetric Drug Discovery Kit

Rapid, sensitive, and simple colorimetric DNMT activity/inhibition kit.

DNMT Colorimetric Drug Discovery Kit is a simple method to detect DNA methylation activity or inhibition with results in <4 hours.

Ordering Information

[Order Online »](#)

ENZ-45017-0001	96Reactions
----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

- Rapid - results in less than 4 hours
- Simple and safe - Easy colorimetric assay without radioactivity or chromatography
- Sensitive - measure as little as 0.5 µg of nuclear extract or 0.5 ng of purified enzymes
- Specific - Optimized to allow high specificity to 5-mC without cross-reactivity to unmethylated cytosine
- Reliable - reproducible results

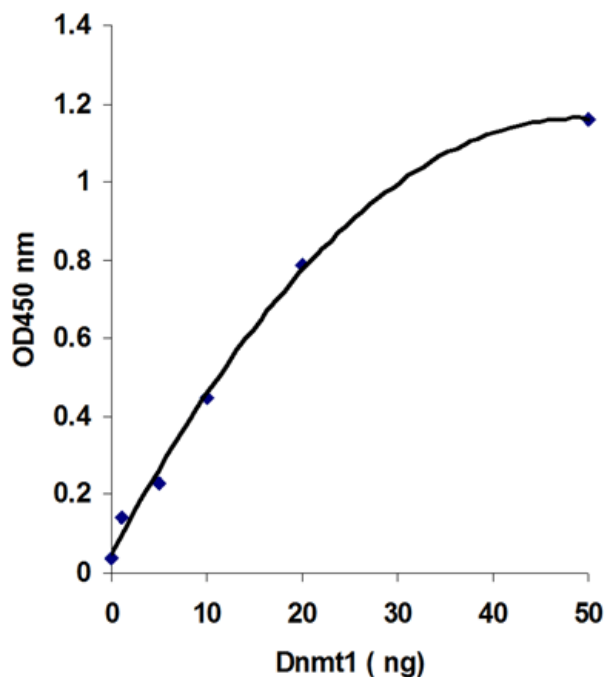


Figure 3. Demonstration of high sensitivity and specificity of DNMT activity/inhibition assay achieved by using recombinant DNMT1 with the DNMT Colorimetric Drug Discovery Kit.

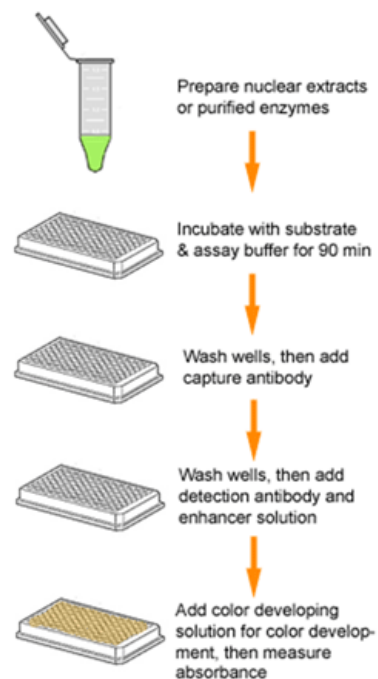


Figure 1. Schematic procedure of the DNMT Colorimetric Drug Discovery Kit.

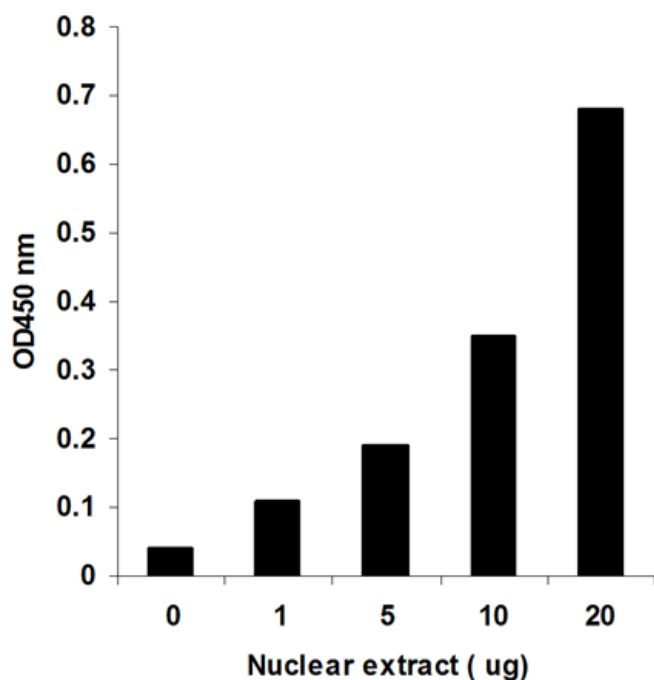


Figure 2. Demonstration of high sensitivity and specificity of the DNMT activity assay achieved by using nuclear extracts with the DNMT Colorimetric Drug Discovery Kit. Nuclear extracts were prepared from MCF-7 cells by using the EPIXTRACT® Nuclear Protein Isolation Kit (Prod. No ENZ-45016).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name DNMT Activity / Inhibition Assay

Application Colorimetric detection

Application Notes For the quantitative determination of DNMT activity/inhibition in nuclear extracts or purified enzymes

Assay Time <4 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Wash Buffer, DNMT Assay Buffer, Adomet, DNMT Enzyme Control, Capture Ab, Detection Ab, Enhancer Solution, Developer Solution, Stop Solution, 8-well Assay Strips with frame, Adhesive Covering Film,



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)