

FLUOR DE LYS[®]- Succinyl Green, Desuccinylase Substrate

The FLUOR DE LYS[®]-*Succinyl Green*, Desuccinylase Substrate offers a highly sensitive and convenient alternative to radiolabeled, succinylated histones or peptide/HPLC methods for the assay of desuccinylases such as SIRT5. The assay procedure has two steps. First, the FLUOR DE LYS[®]-*Succinyl Green*, Desuccinylase Substrate, which comprises a succinylated lysine side chain, is incubated with a sample containing desuccinylase activity (HeLa nuclear or other extract, purified enzyme, bead-bound immunocomplex, etc.). Desuccinylation of the substrate sensitizes the substrate so that, in the second step, treatment with the FLUOR DE LYS[®] Developer (Prod. No. BML-KI105) produces a free rhodamine-based fluorophore which can be monitored with standard fluorometers (Ex. 470-500nm / Em. 520-550nm). The FLUOR DE LYS[®]-*Succinyl Green*, Desuccinylase Substrate has longer wavelength excitation and emission than the AMC-based FLUOR DE LYS[®]-*Succinyl*, Desuccinylase Substrate (Prod. No. BML-KI590). This avoids interference by quenching or fluorescence from compounds absorbing and/or emitting in the near UV and blue range.

Ordering Information

[Order Online »](#)

BML-KI591-0050	50µl
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 1 year when stored at -80°C.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Liquid.
Application Notes	Can be used for the study of desuccinylase kinetics, regulation, and inhibitor sensitivity. Ideal for drug discovery and HTS applications.
Couple Target	HDAC
Couple Type	Substrate
Emission Maximum	520-550nm
Excitation Maximum	470-500nm
Formulation	Solution in DMSO.
Purity	≥95% (TLC)
Source	Synthetic.

Technical Info / Product Notes	NOTE: Must be used in conjunction with FLUOR DE LYS [®] Developer (Prod. No. BML-KI105). For assay conditions, see the product manual for BML-AK514 Green SIRT5 Fluorimetric Drug Discovery Kit.
---------------------------------------	--



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)