## FLUOR DE LYS<sup>®</sup> developer II concentrate

FLUOR DE LYS<sup>®</sup> Developer II is an essential reagent for signal generation in assays done with the FLUOR DE LYS<sup>®</sup>-HDAC8, FLUOR DE LYS<sup>®</sup>-SIRT1 and FLUOR DE LYS<sup>®</sup>-SIRT2 substrates. Deacetylation of the FLUOR DE LYS<sup>®</sup> substrate sensitizes it so that, in a second step, treatment with the FLUOR DE LYS<sup>®</sup> Developer II produces a fluorophore. Typically, an 'inhibitor stop' will be included in the 1x Developer dilution (e.g. nicotinamide for sirtuins).

Citations: 22

View Online »

**Ordering Information** 

Order Online »

**BML-KI176-1250** 1.25ml

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

## **Product Details**

Quantity

NOTE: This product is supplied packaged in 5 tubes (250 µl per tube) for a total volume of 1.25 ml. This packaging is used for optimal stability of the developer.

Last modified: May 29, 2024