

Validation data for Z-IETD-FMK

<https://www.invivogen.com/z-ietd-fmk>

For research use only

Version 21G30-MM

Z-IETD-FMK is a potent small-molecule inhibitor of Caspase-8 (CASP8). Necroptosis, a type of regulated necrosis, is induced when death signals are triggered and CASP-8 is absent or inactivated. The ability of Z-IETD-FMK to induce TNF- α (tumor necrosis factor- α) and RIPK1-dependent necroptosis was validated using InvivoGen's THP1-HMGB1-Lucia™ cellular assay (Figure 1).

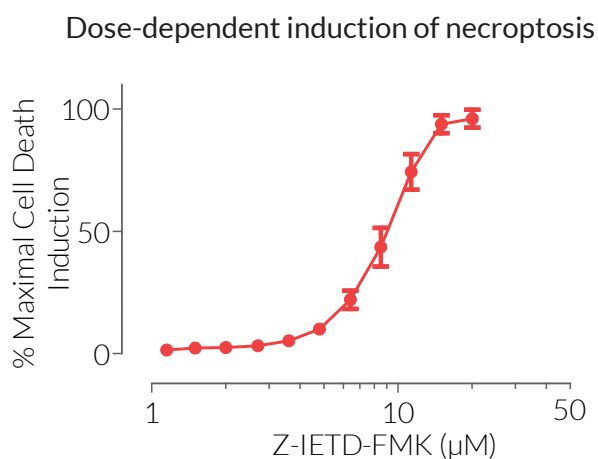


Figure 1: Z-IETD-FMK contributes to TNF- α -induced RIPK1-dependent necroptosis in a dose-dependent manner.

THP1-HMGB1-Lucia™ cells were incubated with recombinant human TNF- α (100 ng/ml), BV6 (cIAP inhibitor, 5 μ M), and increasing concentrations of Z-IETD-FMK (caspase-8 inhibitor). After overnight incubation, the level of HMGB1::Lucia released in the supernatant was assessed by measuring the light signal produced using the QUANTI-Luc™ detection reagent. Data are shown as a percentage (%) of maximal cell death induction.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Asia: +852 3622-3480
E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com