

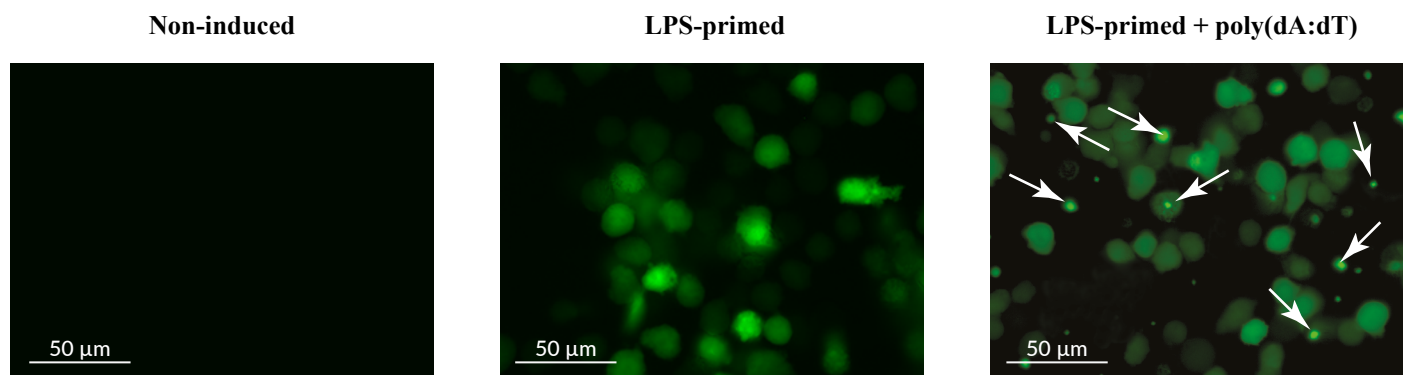
Validation data for THP1-ASC-GFP cells

<http://www.invivogen.com/thp1-asc-gfp>

For research use only

Version 18C14-MM

THP1-ASC-GFP cells are inflammasome reporter cells that allow the monitoring of ASC (apoptosis-associated speck-like protein with a CARD)-dependent inflammasome formation using fluorescence microscopy. These cells stably express a gene encoding an ASC::GFP fusion protein for which expression is driven by an NF- κ B-inducible promoter. Hence, in resting cells, no GFP signal is detected. Upon the first step of inflammasome activation ('priming'), NF- κ B-dependent ASC::GFP expression is induced and can be observed throughout the cytoplasm. Following the second step of inflammasome activation, ASC::GFP polymerizes to form a macromolecular, micrometer-sized complex. The number of ASC::GFP positive cells and localization of fluorescent ASC specks can be determined by using time-lapse confocal or high-resolution fluorescence microscopy.



Visualization of ASC speck formation by fluorescence microscopy: THP1-ASC-GFP cells were primed with 1 μ g/ml LPS-EK for 3 hours, inducing the expression of the ASC::GFP fusion protein in the cytoplasm (middle panel). Cells were then incubated with 250 ng/ml complexed Poly(dA:dT) and ASC speck formation was monitored over 1 to 3 hours post-activation. In most cells, only one speck forms upon inflammasome activation (arrows in right panel). Scale bar: 50 μ m.

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 Hong Kong: +852 3622-3480
E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com