

Validation data for ODN 1826 FITC

<https://www.invivogen.com/odn1826-fitc>

For research use only

Version 24E06-AK

ODN 1826 FITC is a Class B CpG oligonucleotide (ODN). It is a ligand of choice for mouse Toll-like receptor 9 (TLR9). Activation of TLR9 with ODN 1826 triggers NF- κ B- and interferon regulatory factor (IRF)-mediated pro-inflammatory responses. The labeled ligand retains the biological activity of ODN 1826. It can be used to evaluate CpG ODN binding, cellular uptake, and localization by confocal laser-scanning microscopy (excitation 495 nm, emission 520 nm) [data not shown] or flow cytometry (**Figure 1**).

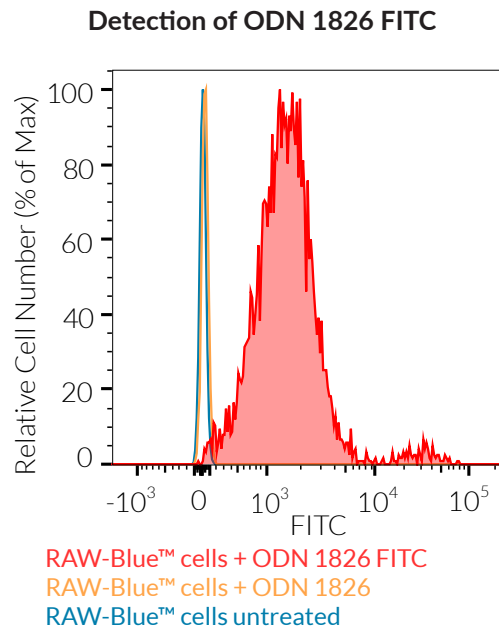


Figure 1. Detection of ODN 1826 FITC. RAW-Blue™ cells were left untreated or incubated with ODN 1826 (unlabeled) or ODN 1826 FITC. After 16h - 24h, the cells were analyzed using flow cytometry.

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