



Focused Screening for Immune-Modulators

- Fully Customizable Compound Screening and Profiling
- Dedicated support team
- Timely and Affordable

Toll-like receptors (TLRs) and other pattern recognition receptors (PRRs) are key targets for drug discovery and vaccine development. With over 25 years of experience in innate immunity, InvivoGen provides tailored compound screening services using highly sensitive, proprietary cell-based assays. Our engineered cell lines are designed to identify lead compounds with immune-modulating properties. Our custom screening and profiling service supports your research in areas such as infectious diseases, cancer, and autoimmune disorders.

Choose from our list of receptor panels:

TLRs (2, 3, 4, 5, 7, 8, 9)

NOD1 and NOD2

Dectin-1, Dectin-2, and Mincle

STING

RIG-I

MDA5

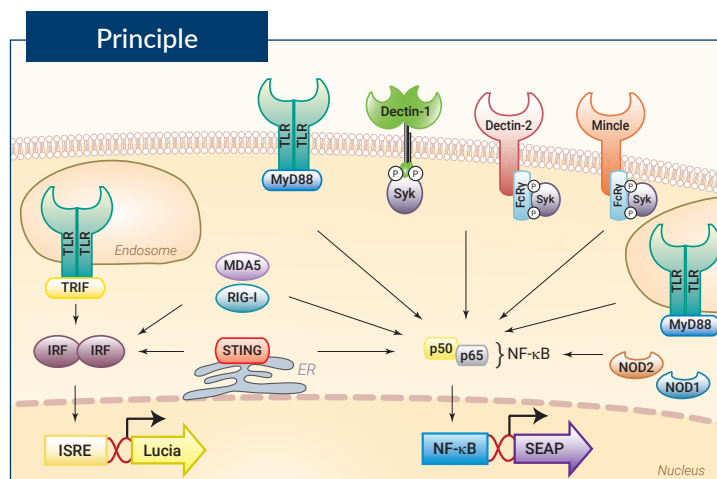
IMMUNOMODULATORY COMPOUND PROFILING SERVICE

InvivoGen's customizable PRR screening services use engineered HEK293, THP-1, or A549 reporter cells to detect activators or inhibitors of key immune pathways. Featuring SEAP and/or Lucia® luciferase reporters to assess NF-κB and IRF pathway activity, our tailored assays deliver precise data, accelerating the discovery of lead compounds.

Our sensitive reporter cell assays offer reliable detection of activators or inhibitors across a comprehensive range of human and mouse pattern recognition receptors:

- **Toll-like Receptors** TLR2, TLR3, TLR4, TLR5, TLR7, TLR8, TLR9
- **NOD-like Receptors** NOD1, NOD2
- **C-type Lectin Receptors** Dectin-1a/b, Dectin-2, Mincle
- **Stimulator of interferon genes (STING)**
- **RIG-I-like Receptors** RIG-I, MDA5

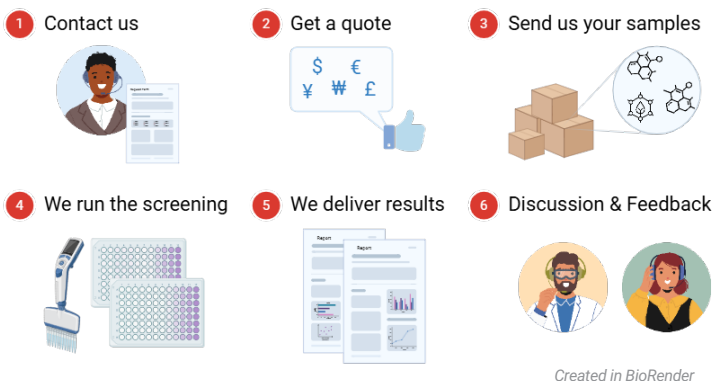
These assays can be performed on all the cell lines in our catalog (e.g. cytokine receptor cells) to meet your specific research needs. Our team will assist you in selecting the most relevant panel for your therapeutic candidates.



Key features

- Fully customizable profiling panels
- Affordable pricing for maximal services
- Suitable for agonist and antagonist screenings
- Fast result delivery (~4 weeks)

Screening service step by step



PRODUCT	CAT. CODE
Compound Profiling	tlrl-test1
Compound Dose Response	tlrl-test2

A detailed report is prepared and provided to the customer electronically. All procedures are performed according to strict guidelines. Confidentiality is guaranteed.

Screening service options

Two choices of services are offered, Compound Profiling (Level 1) and Compound Dose Response (Level 2):

Level 1 - Single dose testing on a set of PRRs

Single-dose testing across one or more selected panels of PRRs to evaluate initial activity. Testing is performed in triplicate.

Level 2 - Dose response on one or several PRRs

Multi-dose testing on specifically chosen PRRs to assess dose dependency. Testing is performed in triplicate.



They trust InvivoGen

More than 60 peer-reviewed publications

Michael J. Newman, Ph.D. Founder and Chief Scientific Officer. Indaptus Therapeutics, Inc.

"I can say without hesitation that InvivoGen is the "go to" source for reagents and research services related to innate immune receptors."

Jeff Meisner, Chief Scientific Officer at Resilient Biotics.

"We've been using InvivoGen products for several years, and their assays consistently deliver high-quality, reproducible data that we can trust."



www.invivogen.com/custom-trl-screening

