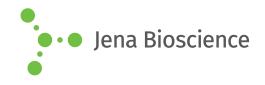
DATA SHEET





NBD-FPP

3,7,11-trimethyl-12-(7-nitro-benzo[1,2,5]oxadiazo-4-ylamino), -dodeca-2,6,10-trien-1 pyrophosphate Fluorescent lipid donor for protein geranylgeranyltransferases (RabGGTase, GGTase-I)

Cat. No.	Amount
LI-013	20 μl

Structural formula of NBD-FPP

For general laboratory use.

Shipping: shipped on gel packs **Storage Conditions:** store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles, store dark

Shelf Life: 12 months

Molecular Formula: C₂₁H₃₀N₄O₁₀P₂ Molecular Weight: 560.43 g/mol

Purity: ≥ 95 % (HPLC)

Form: liquid (Supplied as 1 mM solution in 25 mM (NH₄)₂CO₃)

 $\textbf{Concentration:} \ 1 \ \text{mM}$

Spectroscopic Properties: λ_{exc} 480 nm, λ_{em} 530 nm,

ε 11.0 L mmol⁻¹ cm⁻¹ (pH 7.5)

Description:

NBD-FPP is a fluorescent analog of farnesyl pyrophosphate (FPP). It serves as lipid donor for geranylgeranyltransferases (RabGGTase, GGTase-I). NBD-FPP changes its fluorescence upon binding to protein substrates and allows for efficient fluorescence based GGTase activity assays and inhibitor screening.

Selected References:

Nguyen et al. (2009) Analysis of the eukaryotic prenylome by isoprenoid affinity tagging. Nature Chemical Biology. **5 (4)**:227.

Wu et al. (2007) Synthesis of a fluorescent analogue of geranylgeranyl pyrophosphate and its use in a high-throughput fluorometric assay for Rab geranylgeranyltransferase. Nat. Protoc. 2 (11):2704.

Dursina *et al.* (2006) Identification and specificity profiling of protein prenyltransferase inhibitors using new fluorescent phosphoisoprenoids. *J. Am. Chem. Soc.* **128 (9)**:2822.