

PRODUCT INFORMATION



EYFP mRNA (Cap-1; m1Ψ)

Item No. 43679

Synonym: Enhanced Yellow Fluorescent Protein mRNA (Cap-1; N¹-Methylpseudouridine)

Ex./Em. MAX: 514/527 nm

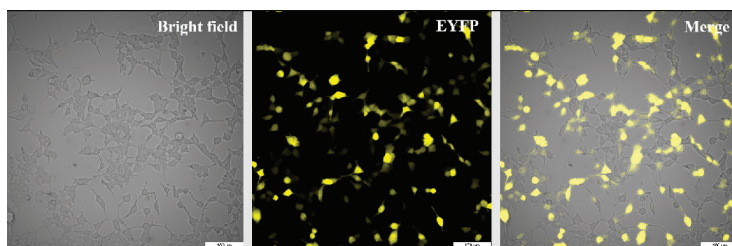
Supplied as: 1 mM sodium citrate, pH 6.4

Storage: -80°C

Stability: ≥6 months

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



EYFP mRNA in 293T cells 24 hours post-transfection.

Description

EYFP mRNA (Cap-1; m1Ψ) encodes for enhanced YFP (EYFP), a yellow fluorescent protein and a derivative of GFP that displays excitation and emission maxima of 514 and 527 nm, respectively.¹ It is capped using a co-transcriptional capping method, resulting in the naturally occurring Cap-1 structure with high capping efficiency. EYFP mRNA is also polyadenylated and modified with N¹-methylpseudouridine (m1Ψ) to reduce host cell immune response and enhance mRNA stability. Encapsulation of EYFP mRNA (Cap-1; m1Ψ) in lipid nanoparticles (LNPs) can be used for mRNA delivery and expression of EYFP.

Reference

1. Suzuki, T., Matsuzaki, T., Hagiwara, H., *et al.* Recent advances in fluorescent labeling techniques for fluorescence microscopy. *Acta Histochem. Cytochem.* **40(5)**, 131-137 (2007).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/05/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM