

# PRODUCT INFORMATION



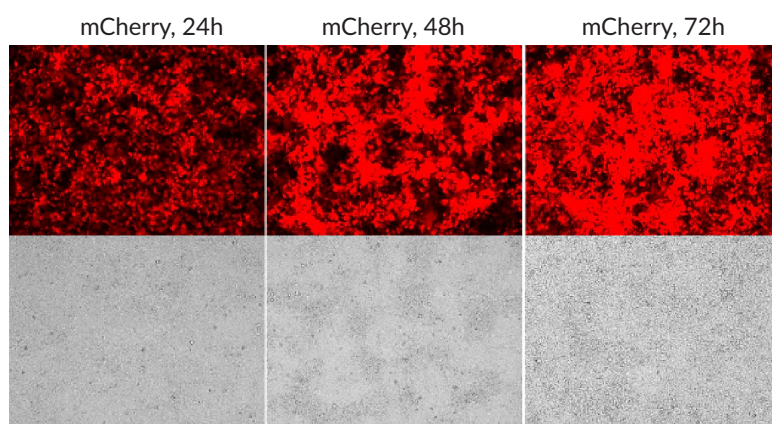
## mCherry circRNA

Item No. 45904

### Overview and Properties

**Synonym:** mCherry Circular RNA  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Supplied as:** A liquid  
**Concentration:** 1 mg/ml  
**Ex./Em. Max:** 587/610 nm

### Image



500 ng circRNA-mCherry was transfected into  $2 \times 10^5$  HEK293 cells by Lip3000 and the brightness of mCherry increased with time after 24h, 48h, and 72h by fluorescence microscopy.

### Description

mCherry circular RNA (circRNA) encodes mCherry, a red fluorescent protein and derivative of *Discosoma* red fluorescent protein (DsRed) that displays excitation/emission maxima of 587/610 nm, respectively.<sup>1</sup> Exogenous circRNAs have increased stability and can produce sustained protein expression compared to mRNA.<sup>2</sup> mCherry circRNA can be used for gene expression, cell labeling *in vitro*, and imaging in animal models.

### References

1. Fernie, A.R. and Willmitzer, L. Molecular and biochemical triggers of potato tuber development. *Plant Physiol.* **127**(4), 1459-1465 (2001).
2. Wesselhoeft, R.A., Kowalski, P.S., and Anderson, D.G. Engineering circular RNA for potent and stable translation in eukaryotic cells. *Nat. Commun.* **9**(1), 2629 (2018).

**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, 05/26/2026

**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