

# PRODUCT INFORMATION



## EZ Cap™ Cre mRNA (ΨUTP)

Item No. 39802

### Overview and Properties

---

**Storage:** -80°C (as supplied)  
**Stability:** ≥6 months  
**Supplied in:** 1 mM Sodium citrate, pH 6.4  
**Concentration:** 1 mg/ml

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

### Description

---

EZ Cap™ Cre mRNA encodes for Cre, a bacteriophage tyrosine recombinase that catalyzes homologous recombination between two loxP sites in DNA.<sup>1</sup> It is capped using a co-transcriptional capping method, resulting in the naturally occurring Cap 1 structure with high capping efficiency. EZ Cap™ Cre mRNA is also polyadenylated and modified with ΨUTP to reduce the host cell immune response and enhance mRNA stability. Encapsulation of EZ Cap™ Cre mRNA (ΨUTP) in lipid nanoparticles (LNPs) can be used for mRNA delivery and expression of Cre *in vitro* or *in vivo*.<sup>2</sup>

### References

---

1. Rajewsky, K., Gu, H., Kühn, R., *et al.* Conditional gene targeting. *J. Clin. Invest.* **98**(3), 600-603 (1996).
2. Kauffman, K.J., Oberli, M.A., Dorkin, J.R., *et al.* Rapid, single-cell analysis and discovery of vectored mRNA transfection *in vivo* with a loxP-flanked tdTomato reporter mouse. *Mol. Ther. Nucleic Acids* **10**, 55-63 (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, 10/12/2023

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