

PRODUCT INFORMATION



LPS from *Salmonella minnesota* R595 (Re)

Item No. 23608

Overview and Properties

Contents: Each vial contains 1 mg of LPS isolated from *Salmonella minnesota* R595 (Re), which was in water when lyophilized

Storage: 4°C (as supplied)

Stability: ≥4 years

Special Conditions: This product is pyrogenic. Avoid accidental autoinoculation by exercising extreme care when handling in conjunction with any injection device.

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

Description

LPS from *Salmonella minnesota* R595 (Re) is a rough mutant strain form of lipopolysaccharide (LPS) extracted from *S. minnesota* R595 (Re). LPS is commonly used to activate toll-like receptor 4 (TLR4) on leukocytes, eliciting inflammatory signaling in isolated cells and *in vivo*.¹⁻⁴

References

1. Mai, K., Chui, J.J., Di Girolamo, N., *et al.* Role of toll-like receptors in human iris pigment epithelial cells and their response to pathogen-associated molecular patterns. *J. Inflamm. (Lond)* **11**, (2014).
2. Medearis, D.N., Camitta, B.M., and Heath, E.C. Cell wall composition and virulence in *Escherichia coli*. *J. Exp. Med.* **128**(3), 399-414 (1968).
3. Migale, R., Herbert, B.R., Lee, Y.S., *et al.* Specific lipopolysaccharide serotypes induce differential maternal and neonatal inflammatory responses in a murine model of preterm labor. *Am. J. Pathol.* **185**(9), 2390-2401 (2015).
4. Pirianov, G., MacIntyre, D.A., Lee, Y., *et al.* Specific inhibition of c-Jun N-terminal kinase delays preterm labour and reduces mortality. *Reproduction* **150**(4), 269-277 (2015).

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, 01/14/2026

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