

PRODUCT INFORMATION



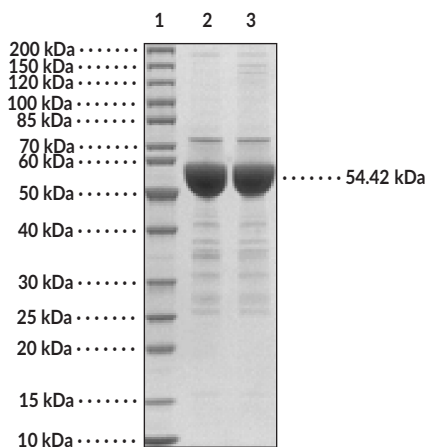
NR6A1 (human, recombinant)

Item No. 41073

Overview and Properties

Synonyms:	GCNF, Germ Cell Nuclear Factor, Nuclear Receptor 61, Nuclear Receptor Subfamily 6 Group A Member 1, Retinoid Receptor-related Testis-associated Receptor
Source:	Recombinant human NR6A1 expressed in <i>E. coli</i>
Amino Acids:	2-480
Uniprot No.:	Q15406
Molecular Weight:	54.42 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Supplied in:	50 mM Tris-HCl, pH 7.5, with 200 mM sodium chloride, 20% glycerol, and 1 mM DTT
Protein Concentration:	<i>batch specific</i> mg/ml

Images



Lane 1: MW Markers
Lane 2: NR6A1 (2 µg, reduced)
Lane 3: NR6A1 (2 µg, non-reduced)

SDS-PAGE Analysis of NR6A1.

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, 07/24/2024

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

PRODUCT INFORMATION



Description

Nuclear receptor subfamily 6 group A member 1 (NR6A1) is a nuclear receptor transcriptional repressor.^{1,2} NR6A1 primarily is expressed in early embryonic development but it is also expressed in developing spermatid cells in testes and pre-meiotic ovaries.³ Nr6a1 homodimers regulate trunk vertebrae number, trunk elongation, and sacral development, as well as promote neuron-related gene expression in neuromesodermal progenitors, during mouse embryonic development.² Deletion of *Nr6a1* is embryonically lethal in mice. Cayman's NR6A1 (human, recombinant) protein consists of 479 amino acids and has a calculated molecular weight of 54.42 kDa.

References

1. Li, J., Mascarinas, P., and McGlenn, E. The expanding roles of Nr6a1 in development and evolution. *Front. Cell Dev. Biol.* **12**, 1357968 (2024).
2. Chang, Y.-C., Manent, J., Schroeder, J., *et al.* Nr6a1 controls *Hox* expression dynamics and is a master regulator of vertebrate trunk development. *Nat. Commun.* **13**(1), 7766 (2022).
3. Chen, F., Cooney, A.J., Wang, Y., *et al.* Cloning of a novel orphan receptor (GCNF) expressed during germ cell development. *Mol. Endocrinol.* **8**(10), 1434-1444 (1994).

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