

PRODUCT INFORMATION



NR4A3 Ligand-binding Domain (human, recombinant)

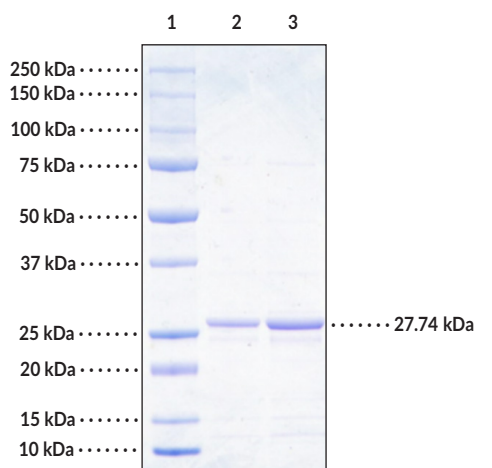
Item No. 40344

Overview and Properties

Synonyms:	Mitogen-induced Nuclear Orphan Receptor, Neuron-derived Orphan Receptor 1, Nuclear Hormone Receptor NOR-1, Nuclear Receptor Subfamily 4 Group A Member 3
Source:	Recombinant human N-terminal His-tagged NR4A3 expressed in <i>E. coli</i>
Amino Acids:	398-626
Uniprot No.:	Q92570
Molecular Weight:	27.74 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥80% estimated by SDS-PAGE
Supplied in:	50 mM Tris HCl, pH 7.5, with 100 mM sodium chloride, and 2% glycerol
Protein Concentration:	<i>batch specific</i> mg/ml

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

Images



Lane 1: MW Markers
Lane 2: NR4A3 Ligand-binding Domain (2 µg)
Lane 3: NR4A3 Ligand-binding Domain (4 µg)

SDS-PAGE Analysis of NR4A3 Ligand-binding Domain.

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, 09/12/2024

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 4 group A member 3 (NR4A3) is a transcription factor and member of the NR4A orphan nuclear receptor family.^{1,2} It is composed of an N-terminal activation function-1 (AF-1) transactivation domain, a DNA-binding domain, a variable linker region, and a C-terminal domain comprising a ligand-binding domain and an AF-2 domain. NR4A3 is primarily expressed in the fetal brain and lungs and adult heart and skeletal muscle, with both splice variants showing similar expression patterns.^{3,4} It is involved in several biological processes, including apoptosis, differentiation, metabolism, inflammation, cardiac remodeling, and development, and its expression can be induced by several factors such as growth factors, cytokines, and stress.^{1,2} Knockout of *Nr4a3* inhibits GM-CSF- and IL-4-induced differentiation of primary mouse monocytes into dendritic cells.⁵ Overexpression of NR4A3 reduces infarct size as a percentage of the area at risk in a mouse model of acute myocardial infarction.⁶ Fusion proteins of NR4A3 with Ewing sarcoma protein (EWS) or TAF15 are associated with extraskeletal myxoid chondrosarcoma.⁷

References

1. Martínez-González, J., Cañes, L., Alonso, J., *et al.* NR4A3: A key nuclear receptor in vascular biology, cardiovascular remodeling, and beyond. *Int. J. Mol. Sci.* **22**(21), 11371 (2021).
2. He, S., Jiang, W., Jiang, B., *et al.* Potential roles of Nr4a3-mediated inflammation in immunological and neurological diseases. *Mol. Neurobiol.* **61**(8), 5958-5973 (2024).
3. Ohkura, N., Ito, M., Tsukada, T., *et al.* Structure, mapping and expression of a human NOR-1 gene, the third member of the Nur77/NGFI-B family. *Biochim. Biophys. Acta* **308**(3), 205-214 (1996).
4. Ohkura, N., Ito, M., Tsukada, T., *et al.* Alternative splicing generates isoforms of human neuron-derived orphan receptor-1 (NOR-1) mRNA. *Gene* **11**(1), 79-85 (1998).
5. Boulet, S., Daudelin, J.-F., Odagiu, L., *et al.* The orphan nuclear receptor NR4A3 controls the differentiation of monocyte-derived dendritic cells following microbial stimulation. *Proc. Natl. Acad. Sci. USA* **116**(30), 15150-15159 (2019).
6. Jiang, Y., Feng, Y.-P., Tang, L.-X., *et al.* The protective role of NR4A3 in acute myocardial infarction by suppressing inflammatory responses via JAK2-STAT3/NF-κB pathway. *Biochem. Biophys. Res. Commun.* **517**(4), 697-702 (2019).
7. Hisaoka, M. and Hashimoto, H. Extraskeletal myxoid chondrosarcoma: updated clinicopathological and molecular genetic characteristics. *Pathol. Int.* **55**(8), 453-463 (2005).

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