

# PRODUCT INFORMATION



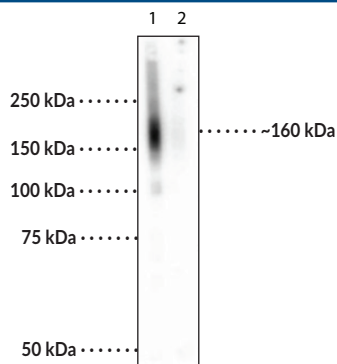
## NCC (Phospho-Thr<sup>53</sup>) Polyclonal Antibody

Item No. 29285

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µl of affinity-purified rabbit polyclonal antibody.
<b>Synonym:</b>	Sodium Chloride Symporter
<b>Immunogen:</b>	Phosphopeptide corresponding to amino acid residues surrounding the phospho-Thr <sup>53</sup> of mouse NCC
<b>Molecular Weight:</b>	~160 kDa
<b>Cross Reactivity:</b>	(+) NCC (phospho-Thr <sup>53</sup> ); (-) Unphosphorylated NCC
<b>Species Reactivity:</b>	(+) Human, mouse, rat
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 µg BSA per ml, and 50% glycerol
<b>Host:</b>	Rabbit
<b>Applications:</b>	Immunofluorescence (IF) and Western blot (WB); the recommended starting dilution is 1:100-1:10,000 and 1:1,000-1:6,000, respectively. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



**Lane 1:** NCC protein phosphorylated at Thr<sup>53</sup>  
**Lane 2:** Phosphospecificity

WB of mouse kidney lysate showing specific immunolabeling of the ~160 kDa NCC protein phosphorylated at Thr<sup>53</sup> in the first lane. Phosphospecificity is shown in the second lane where immunolabeling is completely eliminated by blot treatment with lambda phosphatase (λ-Ptase, 1,200 units for 30 min).

**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, 10/21/2020

**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

---

The thiazide-sensitive sodium chloride cotransporter (NCC) is a member of the SLC12 family of transporters and is encoded by *SLC12A3* in humans.<sup>1</sup> NCC is expressed in epithelial cells of the distal convoluted tubule and localizes to the apical plasma membrane where it mediates sodium reabsorption in the kidney. It consists of a central hydrophobic domain with 12 transmembrane helices that contain affinity-modifying residues for sodium, chloride, and thiazides that is flanked by intracellular N- and C-terminal domains with sites that are subject to phosphorylation. Phosphorylation of NCC at threonine 53 (Thr<sup>53</sup>) is mediated by serum- and glucocorticoid-inducible kinase 1 (SGK1), STE20/SPS1-related proline-alanine-rich protein kinase (SPAK), or oxidative stress-responsive kinase 1 (OSR1). *X. laevis* oocytes expressing a threonine-to-alanine substitution at Thr<sup>53</sup> in NCC, which abolishes its phosphorylation, have reduced chloride deprivation-induced sodium uptake, but not NCC surface expression, compared to wild-type oocytes.<sup>2</sup> NCC (phospho-Thr<sup>53</sup>) levels are increased in the kidney by low dietary intake of sodium chloride or potassium in mice.<sup>3</sup> NCC (phospho-Thr<sup>53</sup>) levels are also increased in the kidney of the CUL3-Het/ $\Delta$ 9 mouse model of familial hyperkalemic hypertension.<sup>4</sup> Cayman's NCC (Phospho-Thr<sup>53</sup>) Polyclonal Antibody can be used for immunofluorescence (IF) and Western blot (WB) applications. The antibody recognizes NCC (Phospho-Thr<sup>53</sup>) at approximately 160 kDa from human, mouse, and rat samples.

## References

---

1. Moes, A.D., van der Lubbe, N., Zietse, R., *et al.* The sodium chloride cotransporter SLC12A3: New roles in sodium, potassium, and blood pressure regulation. *Pflugers Arch.* **466**(1), 107-118 (2014).
2. Pacheco-Alvarez, D., Cristóbal, P.S., Meade, P., *et al.* The Na<sup>+</sup>:Cl<sup>-</sup> cotransporter is activated and phosphorylated at the amino-terminal domain upon intracellular chloride depletion. *J. Biol. Chem.* **281**(39), 28755-28763 (2006).
3. Vallon, V., Schroth, J., Lang, F., *et al.* Expression and phosphorylation of the Na<sup>+</sup>-Cl<sup>-</sup> cotransporter NCC in vivo is regulated by dietary salt, potassium, and SGK1. *Am. J. Physiol. Renal Physiol.* **297**(3), F704-F712 (2009).
4. Ferdaus, M.Z., Miller, L.N., Agbor, L.N., *et al.* Mutant cullin 3 causes familial hyperkalemic hypertension via dominant effects. *JCI Insight* **2**(24), e96700 (2017).

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