

# PRODUCT INFORMATION



## Transcription Factor HIF-1 $\alpha$ Competitor dsDNA

Item No. 10009270

### Overview and Properties

---

<b>Contents:</b>	This vial contains 120 $\mu$ l (10X) specific competitor double-stranded (dsDNA) oligonucleotides containing the HRE sequence (5'-ACGTG-3').
<b>Storage:</b>	-20°C
<b>Stability:</b>	$\geq$ 2 years
<b>Application:</b>	The provided competitor double-stranded oligonucleotide is a specific competitor for HIF-1 $\alpha$ transcription factor binding and demonstrates specificity of the assay. This competitor dsDNA competes with the HRE sequence on the plate for HIF-1 $\alpha$ binding. The competitor oligonucleotide is not biotinylated and therefore will not bind to the plate. In the control nuclear extracts where HIF-1 $\alpha$ specific competitor oligonucleotide has been added, the signal intensity is significantly reduced. The competitor does not need to be run with every assay.

### Procedures

---

Add 10  $\mu$ l of the HIF-1 $\alpha$  specific competitor dsDNA to each well as specified in the HIF-1 $\alpha$  Transcription Factor Assay Kit booklet (Item No. 10006910). These wells should show reduced signal intensity when compared to positive control or sample wells where competitor dsDNA has not been added.

### CAUTIONS

---

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with the HIF-1 $\alpha$  Transcription Factor Assay Kit from Cayman (Item No. 10006910). This product may not perform as described if used with other assay reagents or protocols.

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, 10/18/2021

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