

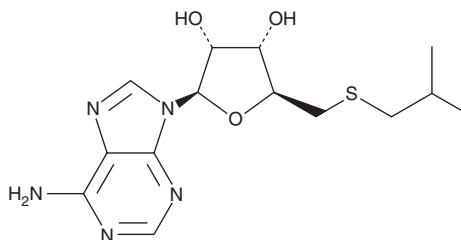
# PRODUCT INFORMATION



## SIBA

Item No. 38758

**CAS Registry No.:** 35899-54-8  
**Formal Name:** 5'-S-(2-methylpropyl)-5'-thio-adenosine  
**Synonyms:** 5'-Deoxy-5'-S-isobutyladenosine, 5'-S-Isobutylthioadenosine  
**MF:** C<sub>14</sub>H<sub>21</sub>N<sub>5</sub>O<sub>3</sub>S  
**FW:** 339.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 261 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

SIBA is supplied as a solid. A stock solution may be made by dissolving the SIBA in the solvent of choice, which should be purged with an inert gas. SIBA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SIBA in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

### Description

SIBA is an isobutylthio-modified derivative of the amino acid derivative S-adenosylhomocysteine (SAH; Item No. 13603).<sup>1</sup> It reduces viral infectivity in HEp-2 cells infected with herpes simplex virus 1 (HSV-1) when used at a concentration of 1 mM.<sup>2</sup> SIBA (0.5 and 1 mM) inhibits the oncogenic transformation of primary chicken embryo fibroblasts induced by Rous sarcoma virus (RSV).<sup>1</sup> It decreases the number of metastases in a model of rhabdomyosarcoma in rats fed a low methionine diet when administered at a dose of 15 mg/kg three times weekly.<sup>3</sup>

### References

1. Robert-Géro, M., Lawrence, F., Farrugia, G., *et al.* Inhibition of virus-induced cell transformation by synthetic analogues of S-adenosyl homocysteine. *Biochem. Biophys. Res. Commun.* **65**(4), 1242-1249 (1975).
2. Jacquemont, B. and Huppert, J. Inhibition of viral RNA methylation in herpes simplex virus type 1-infected cells by 5' S-isobutyl-adenosine. *J. Virol.* **22**(1), 160-167 (1977).
3. Breillout, F., Poupon, M.F., Blanchard, P., *et al.* Association of SIBA treatment and a Met-depleted diet inhibits *in vitro* growth and *in vivo* metastatic spread of experimental tumor cell lines. *Clin. Exp. Metastasis* **6**(1), 3-16 (1988).

#### 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, 07/21/2023

#### 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](http://WWW.CAYMANCHEM.COM)