



## AdenosinePLASMID TAG-LITE® pHALO-A2a Receptor: Adenosine receptor A2a

Product Information Revision: #02 of September 2023

**Part#:** PHTA2A

**Storage temperature:** -20°C

**Quantity:** 10 µg/ 10 µL H<sub>2</sub>O

**Plasmid name:** pHALO-A2a

**Plasmid size:** 7798 bp

**A2A-ORF size:** 2256 bp

**ORF sequence\*:** NM\_000675.4

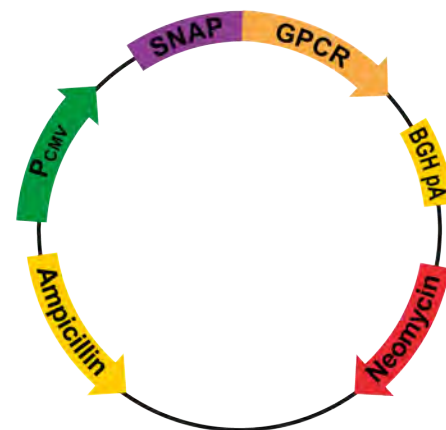
\*based on NCBI nucleic sequence accession number

**Promoter:** CMV

**Host:** Mammalian

**Antibiotic resistance for E. Coli:** Ampicillin

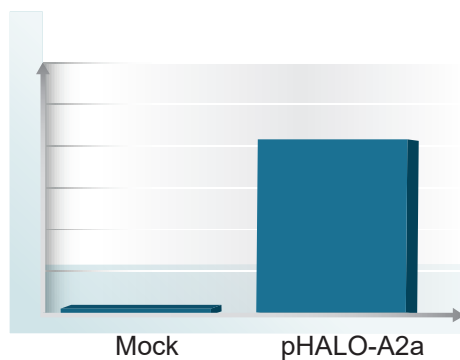
**Mammalian selection marker:** Neomycin



Tag-lite pHALO-A2a

For research use only. Not for use in diagnostic procedures.

### QUALITY CONTROL



Plasmid	Normalized 620 nm signal*
pHALO-A2a	4

\*The plasmid is transiently expressed into HEK 293.

Cells are plated and labelled 24 hours after transfection with the Tag-lite SNAP-Lumi4™-Tb (SSNPTBC) according to the manual described on the web site.

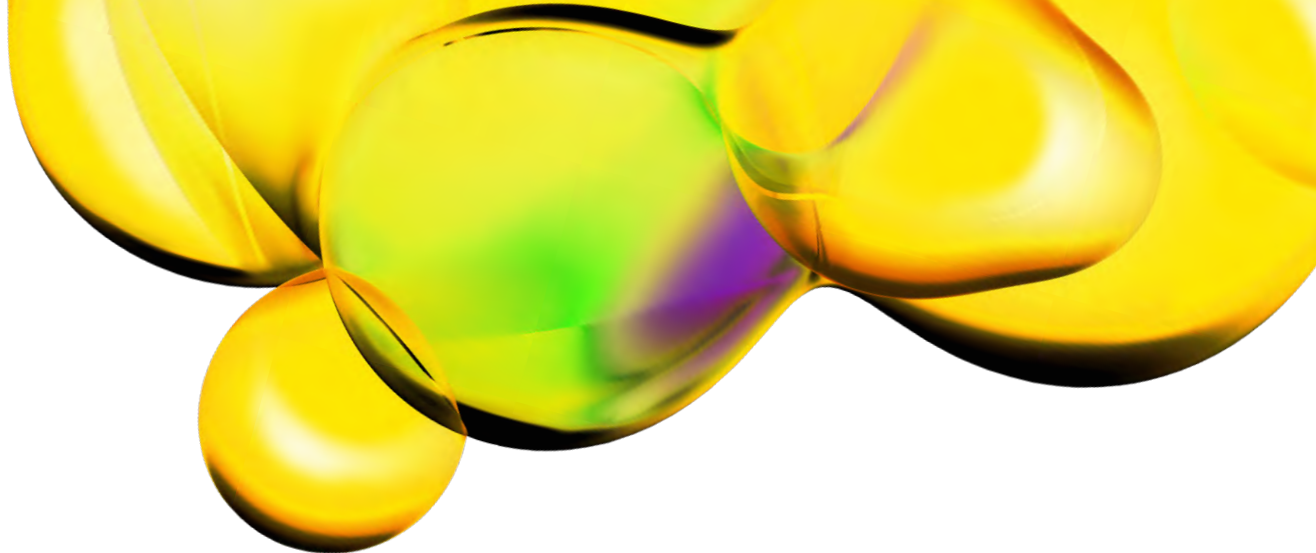
Signal is measured at 620 nm and the value normalized versus the value obtained with a mock plasmid expressed as negative control.

## SEQUENCE & MSDS

MSDS can be downloaded from Tag-lite section on the website [www.revvy.com](http://www.revvy.com).

**Sequence:** M-Halo-H-A2A (923bp - 3178bp, direct) 2256bp

```
ATGGCCTTACCAGTGACCGCCTTGCTCCTGCCGCTGGCCTTGCTGTCTCCACGCCGCCAGGCCGGCCGCCGCTAGCGGCAT
CGAGCAGAAGCTGATCAGCGAGGAGGACCTGGCCGGCATCGATGCAGAAATCGGTACTGGCTTTCCATTTCGACCCCATT
ATGTGGAAAGTCTCGGGCGAGCGCATGCACTACGTCGATGTTGGTCCGCGCGATGGCACCCCTGTGCTGTTCTGCACGGT
AACCCGACCTCCTCCTACGTGTGGCGCAACATCATCCCGCATGTTGCACCGACCCATCGCTGCATTGCTCCAGACCTGAT
CGGTATGGGCAAATCCGACAAACCAGACCTGGGTATTTCTTCGACGACCACGTCCGCTTCATGGATGCCTTCATCGAAG
CCCTGGGTCTGGAAGAGGTGCTCCTGGTCATTACGACTGGGGCTCCGCTCTGGGTTTCCACTGGGCCAAGCGCAATCCA
GAGCGCGTCAAAGGTATTGCATTTATGGAGTTTCATCCGCCCTATCCCGACCTGGGACGAATGGCCAGAATTTGCCCGCA
GACCTTCCAGGCCTTCCGCACCACCGACGTCCGCCGCAAGCTGATCATCGATCAGAACGTTTTTATCGAGGGTACGCTGC
CGATGGGTGTCGTCCGCCCGCTGACTGAAAGTCGAGATGGACCATTACCGCGAGCCGTTCTGAATCCTGTTGACCGCGAG
CCACTGTGGCGCTTCCCAAACGAGCTGCCAATCGCCGGTGAGCCAGCGAACATCGTCGCGCTGGTCCGAAGAATACATGGA
CTGGCTGCACCAGTCCCCTGTCCCGAAGCTGCTGTTCTGGGGCACCCAGGCGTTCTGATCCCACCGGCCGAAGCCGCTC
GCCTGGCCAAAAGCCTGCCTAACTGCAAGGCTGTGGACATCGGCCCGGGTCTGAATCTGCTGCAAGAAGACAACCCGGAC
CTGATCGGCAGCGAGATCGCGCGCTGGCTGTGACGCTGGAGATTTCCGGCGAATTCGATATCCCATCATGGGCTCCTC
GGTGTACATCACGGTGGAGCTGGCCATTGCTGTGCTGGCCATCCTGGGCAATGTGCTGGTGTGCTGGGCCGTGTGGCTCA
ACAGCAACCTGCAGAACGTCACCAACTACTTTGTGGTGTCACTGGCGGGGCCGACATCGCAGTGGGTGTGCTCGCCATC
CCTTTGGCCATCACCATCAGCACCGGGTTCTGCGCTGCCTGCCACGGCTGCCTTTCATTGCCTGCTTCGCTCCTGGTCCT
CACGCAGAGCTCCATCTTCAGTCTCCTGGCCATCGCCATTGACCGCTACATTGCCATCCGCATCCCGCTCCGGTACAATG
GCTTGGTGACCGGCACGAGGGCTAAGGGCATCATTGCCATCTGCTGGGTGCTGTGTTTTGCCATCGGCCTGACTCCCATG
CTAGTTTGAACAACCTGCGGTACGCCAAAGGAGGGCAAGAACCCTCCAGGGCTGCGGGGAGGGCCAAAGTGGCCTGTCT
CTTTGAGGATGTGGTCCCATGAAC TACATGGTGTACTTCAACTTCTTTGCTGTGTGCTGGTGCCCTGCTGCTCATGC
TGGGTGTCTATTTGCGGATCTTCTGGCGGCGGACGACAGCTGAAGCAGATGGAGAGCCAGCCTCTGCCGGGGAGCGG
GCACGGTCCACACTGCAGAAGGAGGTCCATGCTGCCAAGTCACTGGCCATCATTGTGGGGCTCTTTGCCCTCTGCTGGCT
GCCCTACACATCATCAACTGCTTCACTTTCTTCTGCCCCGACTGCAGCCACGCCCTCTCTGGCTCATGTACCTGGCCA
TCGTCCTCTCCACACCAATTCGGTTGTGAATCCCTTCACTACGCCTACCGTATCCGCGAGTTCGCCAGACCTTCCGC
AAGATCATTTCGAGCCACGTCTGAGGCAGCAAGAACCTTTCAAGGCAGTGGCACCAGTGGCCGGGTCTTGGCAGCTCA
TGGCAGTGACGGAGAGCAGGTGAGCCTCCGTCTCAACGGCCACCCGCCAGGAGTGTGGGCCAACGGCAGTGTCCCCACC
CTGAGCGGAGGCCAATGGCTATGCCCTGGGGCTGGTGTGAGTGGAGGGAGTGCCCAAGAGTCCCAGGGGAACACGGGCCTC
CCAGACGTGGAGCTCCTTAGCCATGAGCTCAAGGGAGTGTGCCAGAGCCCCCTGGCCTAGATGACCCCTGGCCCAGGA
TGGAGCAGGAGTGTCC
```



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

[www.revvity.com](http://www.revvity.com)

revvity

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.