

# Substance P (5'- TAMRA conjugate)

Fluorescent photostable ligand for binding to the neurokinin-1 receptor.  
Useful for localizing the NK1R in live tissue slice preparations.

## Ordering Information

[Order Online »](#)

<b>BML-SX5800-0001</b>	1mg
------------------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	TMR-SP, Detection probe for neurokinin-1 receptor
<b>Appearance</b>	Purple powder.
<b>Formulation</b>	Lyophilized.
<b>Identity</b>	Determined by MS.
<b>MW</b>	1761.1
<b>Purity</b>	≥95% (HPLC)
<b>Sequence</b>	Arg-Pro-Lys(5'-TAMRA)-Pro-Gln-Gln-Phe-Phe-Gly-Leu-Met
<b>Solubility</b>	Soluble in water or DMSO (at least 10mg/ml).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)