



## Surelight™ 488 carboxylic acid, NHS ester

Product Number	D44-000-100ug
Amount	100 ug
Molecular Formula	$C_{25}H_{15}Li_2N_3O_{13}S_2$
Molecular Weight	643.41
Au/mg	110.3
Store at	-20°C

### Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at -20°C in the dark.

### Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

<u>Specifications</u>	<u>Actual</u>
Absorbance Max (494 +/- 3nm)	496 nm
Emission Max (518 +/- 4nm)	514.5 nm

For research use only, not for diagnostic or therapeutic use.



**Surelight™ 488 carboxylic acid, succinimidyl ester**

Product Number	D44-000
Amount	1 mg
Molecular Formula	$C_{25}H_{15}Li_2N_3O_{13}S_2$
Molecular Weight	643.41
Au/mg	110.3
Store at	-20°C

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at -20°C in the dark.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

<u>Specifications</u>	<u>Actual</u>
Absorbance Max (494 +/- 3nm)	494 nm
Emission Max (518 +/- 4nm)	516.5 nm

For research use only, not for diagnostic or therapeutic use.

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

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

For technical inquiries: [info@columbiabiosciences.com](mailto:info@columbiabiosciences.com)

For sales inquiries: [sales@columbiabiosciences.com](mailto:sales@columbiabiosciences.com)