

2-Aminoacridone *CAS 27918-14-5*

Catalog Number: 604

Unit Size: 25 mg

Product Details

| | |
|--------------------|--|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
|--------------------|--|

| | |
|-----------------|--------------------------|
| Expiration Date | 12 months upon receiving |
|-----------------|--------------------------|

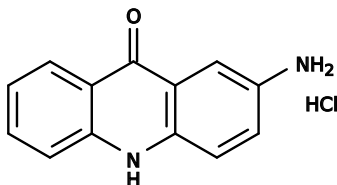
Chemical Properties

| | |
|------------|-------------------|
| Appearance | Solid dark yellow |
|------------|-------------------|

| | |
|------------------|--------|
| Molecular Weight | 246.69 |
|------------------|--------|

| | |
|------------|------|
| Soluble In | DMSO |
|------------|------|

| | |
|--------------------|--|
| Chemical Structure | |
|--------------------|--|

**Spectral Properties**

| | |
|-----------------------|--------|
| Excitation Wavelength | 429 nm |
|-----------------------|--------|

| | |
|---------------------|--------|
| Emission Wavelength | 529 nm |
|---------------------|--------|

Applications

This reagent is used to modify carbonyl-containing biomolecules, in particular, for labeling carbohydrates and glycoproteins. It has large Stokes shift.