

Wheat Germ Agglutinin, XFD488 Labeled

Catalog Number: 25500

Unit Size: 1 mg

Product Details

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

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

Unit Details

| | |
|-------|--------------|
| Units | 25500 (1 mg) |
|-------|--------------|

| | |
|-----------------------|---------------------------|
| Reconstitution Volume | 0.5 mL ddH ₂ O |
|-----------------------|---------------------------|

Chemical Properties

| | |
|------------|--------------|
| Appearance | Solid orange |
|------------|--------------|

| | |
|------------------|-----|
| Molecular Weight | N/A |
|------------------|-----|

| | |
|------------|-------|
| Soluble In | Water |
|------------|-------|

Spectral Properties

| | |
|-----------------------|--------|
| Excitation Wavelength | 499 nm |
|-----------------------|--------|

| | |
|---------------------|--------|
| Emission Wavelength | 520 nm |
|---------------------|--------|

Applications

XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). Wheat germ agglutinin (WGA) is a lectin that binds to N-acetyl-D-glucosamine and sialic acid. It is one of the most studied and useful lectins for its biological applications. Since WGA binds to glycoconjugates its derivatives and conjugates are widely used to label cell membranes and fibrotic scar tissue for fluorescence imaging and analysis. The carbohydrate-binding specificity of WGA is directed against sequences of β -1,4-GlcNAc-linked residues, the chitodextrins. Each monomer contains two identical, non-interacting binding sites, which are complementary to 3 or 4 β -1,4-GlcNAc units. Of the monosaccharides examined, only GlcNAc binds to WGA. ManNAc does not bind, and GalNAc binds only weakly. XFD488 conjugate of WGA is equivalent to the Alexa Fluor® 488 conjugate of WGA. It exhibits bright, green fluorescence. XFD488 WGA conjugate binds to sialic acid and N-acetylglucosaminyl residues.