



Supplier Inclusion Program Frequently Asked Questions

1. Which diverse classifications does Thermo Fisher Scientific accept in the Supplier Inclusion Program?

Our Supplier Inclusion Program provides qualified certified small and diverse businesses with a pathway to become an official provider of goods or services for Thermo Fisher Scientific.

Suppliers that qualify for this program include:

- Minority Owned Business
- Women Owned Business
- LGBT Owned Business
- Veteran Owned Business
- Service-Disabled Veteran Owned Business
- Disabled Owned Business
- Small Businesses registered through the Small Business Administration (SBA)

2. Which certifying agencies does Thermo Fisher Scientific accepts?

Supplier Inclusion Program accepts all State, and local government Supplier Inclusion certifications. In addition, we accept most major national/regional councils such as:

- National Minority Supplier Development Council (NMSDC) www.nmsdc.org
- Women's Business Enterprise National Council (WBENC) www.wbenc.org
- National Veteran-owned Business Association (NaVOBA) www.navoba.com
- National Veteran Business Development Council (NVBDC) www.nvbdc.org
- National Gay and Lesbian Chamber of Commerce (NGLCC) www.nglcc.org
- Disability:IN www.DisabilityIn.org

3. **How does a supplier do business with Thermo Fisher Scientific through its Supplier Inclusion Program?**


To be considered for business through the Supplier Inclusion Program, your company must meet the certification and registration requirements as one of the supplier classifications shown in the previous answer. If your company has met our certification requirements, we encourage you to register with our program through [Gainfront portal](#)



4. What happens after I register with the Supplier Inclusion Program?

After you've completed your registration, your information will reside in our database that is provided to sourcing/procurement staff within Thermo Fisher Scientific. If there's a sourcing opportunity, our buyers and sourcing staff will access this database to search for suppliers based on goods and services, and geographical location information.

5. Who do I contact if I have questions about my STARS (VIVA) registration or technical questions?

- For any technical issues please reach out via email to support@gainfront.com or  [+1-844-717-8815](tel:+18447178815)

6. If I have questions about the Supplier Inclusion Program?

If after reading the Supplier Inclusion Program section of the website you have additional questions please send an email to supplierInclusion@thermofisher.com

