

FAQ for Ductless Air Cleaning Systems

Q: What type of situation would call for the use of the Ductless Air Cleaning System?

A: *When there is an application that requires a high level of filtration, but for various reasons, the application is located in an area without adequate ventilation. The use of the Ductless Air Cleaning System will remove harmful airborne contaminants from the workers breathing zone, and from entering the room environment. Ductless fume hoods are not connected to an exhaust system. They utilize filters to contain vapors, fumes and dusts, before air is recirculated back into the room. Our Ductless Air Cleaning Systems can provide a practical solution for labs where a ducted system is not feasible. Because of their portability, they may be moved from lab to lab and stored away when not in use.*

Q: How do I determine if I need a HEPA Filter Module or a Primary Vapor Removal Module?

A: *Choose the Primary Vapor Removal Modules for chemical fumes and vapors. Choose the Primary HEPA Filter module with vapor after-filter for applications involving powders and particulate such as mold spores and bacteria.*

Q: What type of Filter Media is used in the Primary Vapor Removal Module?

A: *Our Vapor Removal Modules are available with a variety of activated carbons, impregnated carbons, and chemically modified Zeolites. Our special V-Bank design prevents bypass. These filter medias adsorb chemical fumes, including organic solvents, ammonia, acids, and formaldehyde.*

Q: How often should the filter media be replaced?

A: *This is by far the most common question we receive. Every application is different. There are many variables that effect filter life. The type of material being collected, the concentration levels, process temperature, ambient temperature, humidity level, etc. all have an effect on filter life.*

Q: How can I determine when to change the filter media?

A: *Our Ductless Air Cleaning Systems come with a particulate Prefilter. A Prefilter is necessary because it prolongs the life of the HEPA filter. The Prefilter and HEPA filter do not lose efficiency over time. However, they can clog up and reduce the airflow, thus making them less effective over time. When you notice a decrease in performance, first replace the prefilter and then the HEPA filter. The adsorbent media in the Primary Vapor Removal Modules and Vapor Removal After filters should be changed when noticing odor. For most intermittent chemical removal applications, we recommend the filter media be replaced every 3-4 months. If you would like a more specific recommendation please provide details of your specific application to one of our application specialists.*

Q: How Should I dispose of the used filter media?

A: *That depends on the composition of the captured contaminants. Please follow your local regulations.*

Q: What is the motor life expectancy for the Ductless Air Cleaning System?

A: *Our blower motors are totally enclosed and thermally protected. Blower is dynamically balanced. Typical motor life is 100,000 hours, depending on operating conditions, including ambient temperature, duty cycle, etc.*

Q: Can the system be laid on its side to accommodate a height restriction in my space?

A: *Yes, the system can lie horizontally, as well as the typical vertical position. Please make sure to minimize bends or angles in the flexible hose or capture arms when laying the system on its side. Too many bends can lead to restricted airflow.*

Q: There appears to be some loose adsorbent filter media mixed in with the packaging when I first open the box. Is this normal?

A: *Yes, this is normal. During the shipping process, some of the adsorbent media may break down and settle within the module. Because of this, some of the more fine adsorbent material may escape the module. This is completely normal and harmless. Once the module is installed in the system this will no longer occur.*