

Molecular Devices® CellReporterXpress® Software

Third-Party Products Additional Licensing Terms & Conditions

License terms and conditions can be found at the listed website.

Akka.NET	Apache 2 license	https://github.com/akkadotnet/akka.net/blob/dev/LICENSE
Newtonsoft.Json	MIT License	https://github.com/JamesNK/Newtonsoft.Json/blob/master/LICENSE.md
log4net	Apache 2 license	https://svn.apache.org/repos/asf/logging/log4net/tags/log4net-1_2_9/doc/license.html
Castle Core	Apache 2 license	https://github.com/castleproject/Core/blob/master/LICENSE
Nancy	MIT License	https://github.com/NancyFx/Nancy/blob/master/license.txt
Nancy.Authentication.Forms.Owin	MIT License	https://github.com/damianh/Nancy.Authentication.Forms.Owin/blob/master/LICENSE
Nancy.Bootstrappers.Windsor	MIT License	https://github.com/NancyFx/Nancy.Bootstrappers.Windsor/blob/master/license.txt
Nancy.Hosting.Self	MIT License	https://github.com/NancyFx/Nancy/blob/master/license.txt
Nancy.Serialization.JsonNet	MIT License	https://github.com/NancyFx/Nancy.Serialization.JsonNet/blob/master/license.txt
SystemWrapper	Microsoft Public License	https://github.com/jozefizso/SystemWrapper/blob/master/LICENSE.txt
Windsor	Apache 2 license	https://github.com/castleproject/Windsor/blob/master/LICENSE
CsQuery	MIT License	https://github.com/jamietre/CsQuery/blob/master/LICENSE.txt
Iniparser	MIT License	https://github.com/ndevilla/iniparser/blob/master/LICENSE
Topshelf	Apache 2 license	https://github.com/Topshelf/Topshelf/blob/develop/LICENSE
Autofac	MIT License	https://github.com/autofac/Autofac/blob/develop/LICENSE
commonservicelocator	Microsoft Public License	https://github.com/unitycontainer/commonservicelocator/blob/master/LICENSE
Fluent Assertions	Apache 2 license	https://github.com/dennisdoomen/FluentAssertions/blob/master/LICENSE
M2Mqtt	Eclipse Public License 1.0	https://github.com/eclipse/paho.mqtt.m2mqtt/blob/master/LICENSE
NDesk.Options	MIT/X11 license	https://opensource.org/licenses/mit-license.php
Stateless	Apache 2 license	https://github.com/dotnet-state-machine/stateless/blob/dev/LICENSE
OpenSeadragon	New BSD	https://openseadragon.github.io/license/
OpenSeadragonViewerInputHook	MIT License	https://github.com/msalsbery/OpenSeadragonViewerInputHook/blob/master/LICENSE
current-device	MIT License	https://github.com/matthewudson/device.js/blob/master/LICENSE
ag-grid	MIT License	https://github.com/ceolter/ag-grid/blob/master/LICENSE.txt
OpenSeadragonSelection	Apache 2 license	https://github.com/picturae/openseadragonselection/blob/master/LICENSE
angular-busy	MIT License	https://github.com/cgross/angular-busy/blob/master/LICENSE
angular-promise-extras	ISC License	https://github.com/ohjames/angular-promise-extras/blob/master/LICENSE
angular-translate	MIT License	https://github.com/angular-translate/angular-translate/blob/master/LICENSE
angular-tree-control	MIT License	https://github.com/wix/angular-tree-control/blob/master/LICENSE
UI Bootstrap (AngularJS)	MIT License	https://github.com/angular-ui/bootstrap/blob/master/LICENSE
AngularUI Router	MIT License	https://github.com/angular-ui/ui-router/blob/master/LICENSE
animate.css	MIT License	https://github.com/daneden/animate.css/blob/master/LICENSE
bootstrap	MIT License	https://github.com/twbs/bootstrap/blob/v4-dev/LICENSE
interaact.js	MIT License	https://github.com/taye/interact.js/blob/master/LICENSE
jquery	MIT License	http://jquery.org/license
jQuery-Touch-Events	MIT License	https://github.com/benmajor/jQuery-Touch-Events#10-license
ng-onload	MIT License	https://github.com/mikaturunen/ng-onload/blob/master/LICENSE
noUiSlider	MIT License	https://github.com/leongersen/noUiSlider/blob/master/LICENSE.md
ocLazyLoad	MIT License	https://github.com/ocombe/ocLazyLoad/blob/master/LICENSE.md
CsvHelper	Microsoft Public License and Apache 2 license	https://raw.githubusercontent.com/JoshClose/CsvHelper/master/LICENSE.txt
protobuf-csharp-port	(See license file)	https://github.com/jskeet/protobuf-csharp-port/blob/master/license.txt
FileCache	Apache 2 license	https://github.com/acartea/FileCache/blob/master/LICENSE
jose-jwt	MIT License	https://raw.githubusercontent.com/dvsekhvalnov/jose-jwt/master/LICENSE
helios	Apache 2 license	https://github.com/helios-io/helios/blob/master/LICENSE
NLog	BSD 3-Clause "New" or "Revised" License	https://github.com/NLog/NLog/blob/master/LICENSE.txt
SharpZipLib	MIT License	http://www.icsharpcode.net/OpenSource/SharpZipLib/Default.aspx
SQLite.Net-PCL	(See license file)	https://raw.githubusercontent.com/oysteinkrog/SQLite.Net-PCL/master/LICENSE.txt
SQLiteNetExtensions	(See project page)	https://bitbucket.org/twincoders/sqlite-net-extensions
webpack	MIT License	https://github.com/webpack/webpack/blob/master/LICENSE
virtualjoystick.js	MIT License	https://github.com/jeromeetienne/virtualjoystick.js/blob/master/MIT-LICENSE.txt

© 2019 Molecular Devices

V 09/2019