

HSvu Privacy Policy

This Privacy Policy applies to the mobile application HSvu (the "Application") developed by Malvern Panalytical B.V. (the "Service Provider"). The purpose of this policy is to inform users that we do not collect, use, or share any of their personal information.

1. Information Collection and Use

The Application does not obtain or collect any personal or sensitive user information when you download, install, or use it. No registration is required to use the Application, and all data generated during the use of the app is stored exclusively on the user's device and is not transmitted to our servers or any third party. The Application only opens documents stored on your device or shared with you. Furthermore, we do not collect any information on the device, usage data, and location data.

2. Does the Application use third parties?

Since the Application does not collect any information, no data is shared with third parties. The Application does not use analytics tools, advertising networks, external SDKs that collect data.

3. Data Security

Because the Application does not collect or store any personal data, there is no risk of your data being accessed by unauthorized individuals through our service. Data on your device is protected by its operating system's security.

4. Children's Privacy

While the Application is safe for all ages, this Application is not intended for anyone under 13. We do not knowingly collect personal information from children under 13.

5. Changes to This Privacy Policy

We may update this Privacy Policy and will notify you by posting the new policy on this page. Changes are effective immediately upon posting.

This privacy policy is effective as of 14.01.2026.

6. Your Consent

By using the Application, you agree to this Privacy Policy, acknowledging that we do not collect personal information.

7. Contact Us

For questions about this Privacy Policy, contact us at highscore@malvernpanalytical.com