

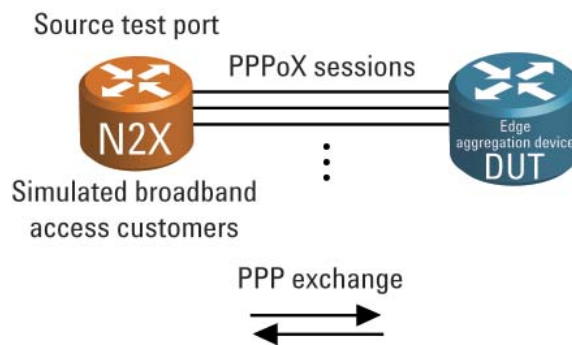
JTC 021 PPPoX Session Setup Rate

Standard RFC 1661, RFC 1332, RFC 1334, RFC 1994.

Purpose This test determines the rate at which a device under test (DUT) can successfully establish “authenticated” PPPoX sessions.

Description In this test, the DUT is an edge aggregation device (e.g., PPPoX server, access concentrator, or router) that supports various access methods, such as PPPoA, PPPoE, and PPPoEoA (referred to generically in this document as “PPPoX”). First, a test port establishes a specified number of PPPoX sessions with the DUT at an initial setup rate. The test then measures the percentage of successful, authenticated sessions the DUT has established, the actual setup time for each session, and an average setup time for that rate. The test is repeated for other setup rates.

Configuration



Procedure

1. From the test port, initiate the establishment of a specified number of PPPoX sessions with the DUT at a specified setup rate (e.g., 25 sessions/second).
2. Measure the number of authenticated sessions the DUT is able to establish and the time it takes to set up each session. Also calculate an average session setup time for that rate.
3. Increase the setup rate on the test port and repeat the test.
4. Continue repeating the test with different setup rates (e.g., 50 sessions/second, 75 sessions/second, 100 sessions/second).

Variables

- Number of PPPoX sessions.
- Setup rates.

Results For each setup rate, record the percentage of successful, authenticated PPPoX sessions the DUT was able to establish as well as the setup times for each session. Also record the average setup time for each setup rate

Graphs Graph the percentage of successful sessions and average setup time for each setup rate, as shown below

